

OptiMAL-IT

Σ 24

REF 710024

Individual Test for Rapid Malaria Diagnostics of *P. falciparum*
et non *P. falciparum* infection in human whole blood



881177 - 2015/03

BIO-RAD

Table of Content

INTRODUCTION	3
REAGENT	3
STORAGE AND VALIDITY OF THE REAGENTS.....	3
WARNINGS AND PRECAUTIONS.....	3
SAMPLE MATERIAL.....	4
POSITIVE CONTROL.....	4
TEST PROCEDURE.....	4
INTERPRETATION OF RESULTS.....	5
REMARK.....	6
LIMITATIONS.....	6
BIBLIOGRAPHY.....	6
PRODUCT.....	6

INTRODUCTION

OptiMAL-IT is an immuno-chromatographic test, using monoclonal antibodies against the metabolic enzyme pLDH (parasite lactate dehydrogenase) of *Plasmodium spp.* These monoclonal antibodies are classified in two groups:

- 1) one specific for *Plasmodium falciparum*
- 2) and the other is a pan-specific monoclonal antibody which reacts with all four species of *Plasmodium spp.* which can occur in human beings: *P. falciparum*, *P. vivax*, *P. ovale*, *P. malariae*.

In case of presence of *Plasmodium spp.* in the blood sample, the pLDH captured by the conjugate reacts with the specific antibodies against *Plasmodium falciparum* and/or *Plasmodium spp.*

Specificity: **OptiMAL-IT** detects the presence of pLDH, an enzyme produced by both sexual and asexual form of the parasite. There is no cross-reaction with human LDH or Rheumatoid factor. The “lysis and dilution” buffer avoids also non-specific reactions due to heterophiles antibodies.

Sensitivity: **OptiMAL-IT** can detect peripheral parasitaemia levels of 0.001–0.002% (50-100 parasites per μL of blood). This sensitivity can be compared to microscopic observation of a thin blood smear using a x100 immersion objective, for a period of approx. 30 minutes, by a well-trained technician.

Fast and easy: The result is obtained in 20 minutes. **OptiMAL-IT** provides all necessary materials to perform a test.

Monitoring of treatment with OptiMAL-IT: the test is only positive when live parasites are present in the blood. A repeat test becomes negative generally within 2-4 days following the beginning of successful treatment. Thus **OptiMAL-IT** is suitable to verify the effectiveness of therapy, underlining possible resistant strains of *Plasmodium spp.*

REAGENT

Each single test package contains:

- 1 device with dipstick, conjugate well, wash well
- 1 well cover
- 1 dropper ampoule with buffer
- 1 lancet
- 1 disinfecting swab
- 1 pipette (printed mark for 10 μL)
- 1 schematic test procedure.

STORAGE AND VALIDITY OF THE REAGENTS

OptiMAL-IT must be stored at +2-30°C. When stored in correct conditions, **OptiMAL-IT** is stable until the expiry date on aluminium foil. Do not freeze.

Important: When transporting or storing the packages, avoid exposure to high temperature (over 40 °C) for a period longer than 1 day. Avoid any exposure to 60 °C and higher (the reagents may be damaged).

WARNINGS AND PRECAUTIONS

Reliability of results depends on correct implementation of Good Laboratory Practices.

- For *in vitro* diagnostic use only.
- Do not use any part of the test beyond the expiration date.
- Always follow recommended storage conditions.
- For single use only.
- Handle patient samples and any material that comes directly in contact with them as potentially able to transmit infectious diseases.

- Wear disposable gloves when handling samples and reagents.
- Do not touch the reaction zone with fingers.
- Avoid spilling samples. Spills must be rinsed with bleach diluted to 10 %. The material used for cleaning must be discarded in a contaminated residue container.
- Patient samples and contaminated material should be discarded after decontamination.
- For hazard and precaution recommendations related to some chemical components in this test kit, please refer to the pictogram(s) mentioned on the labels and the information supplied at the end of instruction for use. The Safety Data Sheet is available on www.bio-rad.com.

SAMPLE MATERIAL

- Capillary blood collected from fingertip.
- Whole blood collected by venipuncture, using EDTA sample tubes.

POSITIVE CONTROL (separately available)

OptiMAL Positive control is a recombinant pLDH giving exactly the same result as a blood sample from a patient infected with *Plasmodium falciparum*.

It is recommended to use this reagent to verify the performance of **OptiMAL-IT** in a regular manner, and especially after transportation and storage under tropical conditions.

For use, consult the specific package insert.

TEST PROCEDURE

Tear open the aluminium package and take out all the material.

Important: *Do not leave the material exposed to humidity and high temperature. In tropical conditions, use the test within 15 minutes after opening the aluminium package.*

1. Take the device, place it horizontally on a flat surface, write the patient's name or number on the label.
2. Tear open the ampoule of buffer, add **1 drop** of buffer to the first well (conjugate well, marked with a colored line), and **4 drops** to the second well (wash well). Allow to stand for **1 minute**.
3. For finger prick blood: clean the fingertip with the swab, let dry, remove the lancet from its envelope, prick the finger. Take the pipette, squeeze it, place the open tip into the blood drop, release pressure and draw up blood to the black line. Discard used swab and lancet into a suitable waste container. When using venous blood, draw blood from the tube into the pipette in the same manner.
4. Add the entire volume of blood by squeezing the pipette gently, to the first well (conjugate well, marked with a colored line).
5. Stir gently with the upper end of the pipette and allow to stand for **1 minute**. Discard the pipette into a suitable waste container.
6. Pull the IT device apart: hold the device with the wells between thumb and forefinger and, with the other hand, pull out the dipstick holder (with the label). Place the wells back on the table, insert the legs of the dipstick holder into the holes beside the conjugate well (with colored line) so that the dipstick end reaches the bottom of the conjugate well. Allow to stand for **10 minutes**. The blood/conjugate mixture should then be completely soaked up.
7. Transfer the dipstick to the second well (wash well) and allow to stand for **10 minutes**. The reaction field should then be completely cleared of blood. The control band must be clearly visible.
8. Remove the dipstick from the wash well and click it back into the clear plastic piece. Close the wells with the well cover, break them off, and break the two legs off from the clear plastic piece. Discard them into a suitable waste container.
9. Read the reaction and interpret the results (see: Interpretation). The dipstick slide should be kept for future reference and for comparison in case of further tests for monitoring the efficiency of treatment.

For routine use, you may follow the schematic test procedure enclosed in each test package.

INTERPRETATION OF RESULTS

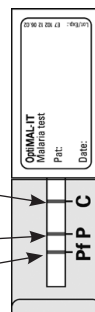
A) Principle

OptiMAL-IT is a rapid detection test using a dipstick coated with monoclonal antibodies against the metabolic enzyme pLDH (parasite lactate dehydrogenase).

In case of presence of *Plasmodium* in the blood sample, the pLDH captured by the conjugate reacts with the specific antibodies against *Plasmodium falciparum* and/or *Plasmodium* spp.

The reactions are demonstrated by the appearance of dark purple bands on the dipstick. The band of the procedure control must always be present to validate a result. In case of absence of the control band, the test is not valid.

- Goat anti-mouse (procedure control) «C»
- Monoclonal anti-pLDH reacts with all 4 species of *Plasmodium*:
P. falciparum / *P. vivax* / *P. malariae* / *P. ovale* «P»
- Monoclonal anti-pLDH specific only to *P. falciparum* «Pf»



B) Validation of the test

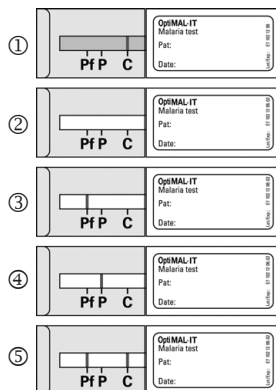
Results are valid if:

- the control band is clearly visible.
- the reaction field is cleared of blood.

Results are NOT valid if:

- the dipstick is not sufficiently cleared (reaction field remains red) [1].
- the control band is not present [2].
- the control band is not visible even if one or both of the diagnostic bands are present [3][4].
- the control band and the band «Pf» (*P. falciparum*) are present while the band «P» is absent [5].

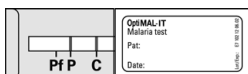
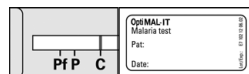
In such cases, repeat the test, following precisely the test procedure.



C) Interpretation of the reaction pattern

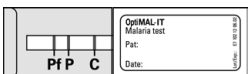
Negative reaction: No detectable pLDH is present in the sample.

Positive reaction: The pLDH present in the sample reacts with the anti-pLDH conjugate and rises up the Dipstick where it is captured by one or both of the specific pLDH antibodies.



POSITIVE for *P. vivax*, or *P. malariae*, or *P. ovale*.

This sample is **negative** for *P. falciparum*.



POSITIVE for *P. falciparum*

Remark: In case of a positive reaction for *Plasmodium falciparum*, the possibility of a mixed infection with *Plasmodium vivax*, *P. ovale* or *P. malariae* cannot be excluded.

D)Monitoring of treatment

The test is only positive when live parasites are present in the blood; thus, **OptiMAL-IT** is also suitable to verify the effectiveness of therapy, underlining possible resistant strains of *Plasmodium spp.*

2 days and again 4-5 days after start of the treatment repeat testing and compare the reactions with the results obtained before treatment.

Weaker positive reactions after 2 days and negative reaction after 4-5 days indicate an efficient treatment.

If the reaction remains positive after 5-7 days, the possibility of a resistant strain should be considered.

Depending on the medication, the treatment may take longer. In this case, repeat the test 7 to 10 days after the start of treatment.

REMARK

The unused dipsticks present faint lines at the position of the 3 antibodies. These faint lines are present for the quality control during the manufacturing process. They disappear completely as soon as the reaction field of the dipstick is in contact with the blood sample and conjugate.

LIMITATIONS

- **OptiMAL-IT** detects the presence of pLDH from live parasites only (hence the possibility to monitor the efficiency of treatment). Comparison tests with PCR may therefore not be conclusive.
- Any modification of the described test procedure or use of other reagents may modify the reaction pattern and invalidate the test.
- Remember that the result of **OptiMAL-IT** is to be interpreted within the epidemiological, clinical and therapeutical context.

BIBLIOGRAPHY

1. Lactate Dehydrogenase and the Diagnosis of Malaria. Makler M.T., Piper R.C., & Milhous. Parasitology Today, 14(9), 376-377, 1998.
2. Performance of the OptiMAL malaria antigen capture dipstick for malaria diagnosis and treatment monitoring at the Hospital for Tropical Diseases, London. Moody A.H., Hunt Cooke A., Gabbett E. and Chodini P. Br.J.Haematol., 109, 891-894, 2000.
3. Malaria diagnosis and treatment; relevance of laboratory support from the rapid immunochromatographic tests of ICT-Malaria (P.f/P.v) and OptiMAL. Tarimo DS., Minjas JN., Bygbjerg IC. Ann.Trop.Med.Parasit., 95[5], 437-444, 2001.
4. Rapid Diagnostic Tests for Malaria Parasites. A. Moody. Clin. Microbiol. Rev. 15 [1], 66-78, 2002.
5. Non-microscopic method for malaria diagnosis using OptiMAL-IT, a second-generation dipstick for malaria pLDH antigen detection. Moody A. H. and Chiodini P. L. British Journal of Biomedical Science, 59 [4], 228-321, 2002.

PRODUCT

OptiMAL-IT Individual Rapid Malaria test Box of 24 testsREF 710024
OptiMAL Positive control 1 Tube of 8 wellsREF 710011

- (BG)** • Този продукт съдържа човешки или животински компоненти. Бъдете внимателни при работа с него.
- (CZ)** • Tento výrobek obsahuje lidské nebo zvířecí komponenty. Zacházejte s ním opatrně.
- (DE)** • Dieses Produkt enthält Bestandteile menschlichen oder tierischen Ursprungs. Vorsichtig handhaben.
- (DK)** • Dette produkt indeholder humane og animalske komponenter. Skal behandles med forsigtighed.
- (EE)** • Käesolev toode sisaldab inim-või loomseid komponente. Käsitseta ettevaatlikult.
- (EN)** • This product contains human or animal components. Handle with care.
- (ES)** • Este producto contiene componentes humanos o animales. Manejar con cuidado.
- (FI)** • Tässä tuotteessa on ihmisestä tai eläimestä peräisin olevia osia. Käsittele varovasti.
- (FR)** • Ce produit contient des composants d'origine humaine ou animale. Manipuler avec précaution.
- (GR)** • Αυτό το προϊόν περιέχει ανθρώπινα ή ζωικά στοιχεία. Χειριστείτε το με προσοχή.
- (HR)** • Ovaj proizvod sadrži ljudske ili životinjske sastojke. Pažljivo rukovati.
- (HU)** • A készítmény emberi vagy állati eredetű összetevőket tartalmaz. Óvatosan kezelendő.
- (IT)** • Questo prodotto contiene componenti umane o animali. Maneggiare con cura.
- (LT)** • Šiame produkto yra žmogiškosios arba gyvūninės kilmės sudėtiniai dalys. Elgtis atsargiai.
- (NL)** • Dit product bevat menselijke of dierlijke bestanddelen. Breekbaar.
- (NO)** • Dette produktet inneholder humane eller animalske komponenter. Håndteres med forsiktighet.
- (PL)** • Niniejszy produkt zawiera składniki pochodzenia ludzkiego lub zwierzęcego. Należy obchodzić się z nim ostrożnie.
- (PT)** • Este medicamento contém componentes de origem humana ou animal. Manuseie com cuidado.
- (RO)** • Acest produs conține materiale de origine umană sau animală. Manevrați-l cu grijă.
- (SE)** • Denna produkt innehåller beståndsdelar från människa eller djur. Hantera produkten varsamt.
- (SI)** • Izdelek vsebuje človeške ali živalske sestavine. Rokujte previdno.
- (SK)** • Tento výrobok obsahuje ľudské alebo zvieracie zložky. Narábajte s ním opatrne.



H360-H319
P280-P337+P313
P305+P351+P338-P501

(CZ)

Varování

Může poškodit reprodukční schopnost nebo plod v těle matky. Způsobuje vážné podráždění očí.
Používejte ochranné rukavice/ochranný oděv/ochranné brýle/obličejový štít. Přežívá-li podráždění očí:
Vyhleďte lékařskou pomoc/ošetření. **PŘI ZASAŽENÍ OČÍ:** Několik minut opatrně vyplachujte vodou. Vyměte kontaktní čočky, jsou-li nasazeny a pokud je lze vyměnit snadno. Pokračujte ve vyplachování. Obsah/nádobu likvidujte v souladu s místními/regionálními/národními/mezinárodními předpisy.

(DE)

Achtung

Kann die Fruchtbarkeit beeinträchtigen oder das Kind im Mutterleib schädigen. Verursacht schwere Augenreizung. Schutzhandschuhe/Schutzkleidung/Augenschutz/ Gesichtsschutz tragen. Bei anhaltender Augenreizung: Ärztlichen Rat einholen/ärztliche Hilfe hinzuziehen. **BEI KONTAKT MIT DEN AUGEN:** Einige Minuten lang behutsam mit Wasser spülen. Vorhandene Kontaktlinsen nach Möglichkeit entfernen. Weiter spülen. Entsorgung des Inhalts / des Behälters gemäß den örtlichen / regionalen / nationalen/ internationalen Vorschriften.

(DK)

Advarsel

Kan skade forplantningsevnen eller det ufødte barn. Forårsager alvorlig øjenirritation. Bær beskyttelsehandsker/beskyttelsestøj/ øjenbeskyttelse/ansigtsbeskyttelse Ved vedvarende øjenirritation: Søg lægehjælp. **VED KONTAKT MED ØJNENE:** Skyl forsigtigt med vand i flere minutter. Fjern eventuelle kontaktlinser, hvis dette kan gøres let. Fortsæt skylning. Bortskaffelse af indholdet/holderen i henhold til de lokale/regionale/nationale/internationale forskrifter.

(EE)

Hoiatus

Võib kahjustada viljakust või loodet. Põhjustab tugevat silmade ärritust.

Kanda kaitsekindaid/kaitserõivastust/kaitseprille/ kaitsemaski. Kui silmade ärritus ei möödu: pööruda arsti poole. **SILMA SATTUMISE KORRAL:** loputada mitme minuti jooksul ettevõtteilikut veega. Eemaldada kontaktläätsed, kui neid kasutatakse ja kui neid on kerge eemaldada. Loputada veel kord. Sisukonteineri käitlus vastavuses kohalike/regionaalsete/rahvuslike/ rahvusvaheliste nõuetega.

(EN)

Warning

May damage fertility or the unborn child. Causes serious eye irritation.

Wear protective gloves/protective clothing/eye protection/ face protection. If eye irritation persists: Get medical advice/attention. **IF IN EYES:** Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Dispose of contents/ container in accordance with local/regional/national/ international regulations.

(ES)

Atención

Puede perjudicar la fertilidad o dañar al feto. Provoca irritación ocular grave. Llevar guantes que aislen del frío/gafas/máscara. Si persiste la irritación ocular: Consultar a un médico. **EN CASO DE CONTACTO CON LOS OJOS:** Aclarar cuidadosamente con agua durante varios minutos. Quitar las lentes de contacto, si lleva y resulta fácil. Seguir aclarando. Eliminar el contenido o el recipiente conforme a la reglamentación local/regional/nacional/ internacional.

(FI)

Varoitus

Saattaa heikentää hedelmällisyyttä tai vaurioittaa sikiötä. Ärsyttää voimakkaasti silmiä. Käytä suojäkäsineitä/suojavaateetusta/silmien suojausta/ kasvonsuojausta. Jos silmää ärsytys jatkuu: Hakeudu lääkäriin. **JOS KEMIKAALIA JOUTUU SILMIIN:** Huuhto huolellisesti vedellä usean minuutin ajan. Poista piilolinssit, _edical voi tehdä helposti. Jatka huuhtomista. Säilytä säiliö(t) noudattaen paikallisia/alueellisia/ kansallisia/kansainvälisiä määräyksiä.

(FR)

Attention

Peut nuire à la fertilité ou au fœtus. Provoque une sévère irritation des yeux. Porter des gants de protection/des vêtements de protection/un équipement de protection des yeux/ du visage. Si l'irritation oculaire persiste: consulter un médecin. **EN CAS DE CONTACT AVEC LES YEUX:** rincer avec précaution à l'eau pendant plusieurs minutes. Enlever les lentilles de contact si la victime en porte et si elles peuvent être facilement enlevées. Continuer à rincer. Éliminer le contenu/réceptacle conformément à la réglementation locale/régionale/nationale/internationale.

(GR)

Προσοχή

Μπορεί να βλάψει τη γονιμότητα ή το έμβρυο. Προκαλεί σοβαρό οφθαλμικό ερεθισμό. Να φοράτε προστατευτικά γάντια/προστατευτικά ενδύματα/μέσα ατομικής προστασίας για ταμάτια/ πρόσωπο. Εάν δεν υποχωρεί ο οφθαλμικός ερεθισμός: Συμβουλευθείτε/Επισκεφθείτε γιατρό. **ΣΕ ΠΕΡΙΠΤΩΣΗ ΕΠΑΦΗΣ ΜΕ ΤΑ ΜΑΤΙΑ:** Ξεπλύνετε προσεκτικά με νερό για αρκετά λεπτά. Εάν υπάρχουν φακοί επαφής, αφαιρέστε τους, εφόσον είναι εύκολο. Συνεχίστε να ξεπλύνετε. Απορρίψτε τα περιεχόμενα/δοχείο σύμφωνα με τους τοπικούς/εθνικούς/διεθνείς κανονισμούς.

(HR)

Upozorenje

Može štetno djelovati na plodnost ili naškoditi nerođenom djetetu. Uzrokuje jako nadraživanje oka. Nositi zaštitne rukavice/zaštitnu odijelo/zaštitu za oči/ zaštitu za lice. Ako nadražaj oka ne prestaje: zatražiti savjet/pomoć liječnika. **U SLUČAJU DODIRA S OČIMA:** oprezno ispirati vodom nekoliko minuta. Ukloniti kontaktne leće ukoliko ih nosite i ako se one lako uklanjaju. Nastaviti ispiranje. Odložite sadržaje / spremnike u skladu s lokalnim/regionalnim/nacionalni/ međunarodnim odredbama.

(HU)

Figyelem

Károsíthatja a termékenységet vagy a születendő gyermeket. Súlyos szemirritációt okoz.

Védőkesztyű/védőruha/szemvédő/arcvédő használata kötelező. Ha a szemirritáció nem múlik el: orvosi ellátást kell kérni. SZEMBE KERÜLÉS esetén: Több percig tartó óvatos öblítés vízzel. Adott esetben a kontaktlencsék eltávolítása, ha könnyen megoldható. Az öblítés folytatása. Az edény tartalmát / a tartályt a helyi/regionális/nemzeti/nemzetközi szabályozásoknak megfelelően kell hulladékként elhelyezni.

(IT)

Attenzione

Può nuocere alla fertilità o al feto. Provoca grave irritazione oculare.

Indossare guanti/indumenti protettivi/Proteggere gli occhi/ il viso. Se l'irritazione degli occhi persiste, consultare un medico. IN CASO DI CONTATTO CON GLI OCCHI: sciacquare accuratamente per parecchi minuti.

Togliere le eventuali lenti a contatto se è agevole farlo. Continuare a sciacquare. Smaltire il prodotto/recipiente in conformità con le disposizioni locali / regionali / nazionali / internazionali.

(LT)

Atsargiai

Gali pakenkti vaisingumui arba negimusiam vaikui.

Sukelia smarkų akių dirginimą.

Mūvėti apsaugines pirštines/dėvėti apsauginius drabužius/naudoti akių (veido) apsaugos priemones. Jei akių dirginimas nepaėina: kreiptis į gydytoją. PATEKUS Į AKIS: Kelias minutes atsargiai plauti vandeniu. Išimti kontaktinius lęšius, jeigu jie yra ir jeigu lengvai galima tai padaryti. Toliau plauti akis. Turinį/talpą išpilti (išmesti) - šalinti pagal vietines / regionines / nacionalines / tarptautines taisykles.

(NL)

Waarschuwing

Kan de vruchtbaarheid of het ongeboren kind schaden. Veroorzaakt ernstige oogirritatie.

Beschermende handschoenen/beschermende kleding/oogbescherming/gelaatsbescherming dragen. Bij aanhoudende oogirritatie: een arts raadplegen. BIJ CONTACT MET DE OGEN: voorzichtig afspoelen met water gedurende een aantal minuten; contactlenzen verwijderen, indien mogelijk; blijven spoelen. De inhoud en de verpakking verwerken volgens de plaatselijke/regionale/nationale/internationale voorschriften.

(NO)

Advarsler

Kan redusere forplantningsevnen eller gi fosterskader.

Forårsaker alvorlige øyeirritasjoner.

Bruk vernehansker/vermekåper/vernebriller/ansiktsskjerm. Ved vedvarende øyeirritasjon: Kontakt / tilkall lege. VED KONTAKT MED ØYNENE: Skyll forsiktig med vann i opptil flere minutter. Fjern evt. kontaktlinser såfremt dette er lett mulig. Fortsett skyllingen. Innholdet / emballasjen skal avhendes i henhold til de lokale / regionale / nasjonale / internasjonale forskrifter.

(PL)

Uwaga

Może działać szkodliwie na płodność lub na dziecko w łonie matki. Działa drażniąco na oczy.

Stosować rękawice ochronne/odzież ochronną/ochronę oczu/ochronę twarzy. W przypadku utrzymywania się działania drażniącego na oczy: Zasięgnąć porady/zgłosić

się pod opiekę lekarza. W PRZYPADKU DOSTANIA SIĘ DO OCZU: Ostrożnie płukać wodą przez kilka minut. Wyjąć soczewki kontaktowe, jeżeli są i można je łatwo usunąć. Nadal płukać. Zawartość / pojemnik usuwać zgodnie z przepisami miejscowymi / regionalnymi / narodowymi / międzynarodowymi.

(PT)

Atenção

Pode afectar a fertilidade ou o nascituro. Provoca irritação ocular grave.

Usar luvas de protecção/vestuário de protecção/protecção ocular/protecção facial. Caso a irritação ocular persista: consulte um médico. SE ENTRAR EM CONTACTO COM OS OLHOS: enxaguar cuidadosamente com água durante vários minutos. Se usar lentes de contacto, retire-as, se tal lhe for possível. Continuar a enxaguar. Eliminar o conteúdo/recipiente de acordo com a legislação local/regional/nacional/internacional.

(RO)

Atenție

Poate dauna fertilității sau fătului. Provoacă o iritare gravă a ochilor.

Purtați mănuși de protecție/îmbrăcămintele de protecție/echipament de protecție a ochilor/ chipament de protecție a feței. Dacă iritarea ochilor persistă: consultați medicul. ÎN CAZ DE CONTACT CU OCHI: clătiți cu atenție cu apă timp de mai multe minute. Scoateți lentilele de contact, dacă este cazul și dacă este lucru se poate face cu ușurință. Continuați să clătiți. Aruncați conținutul/ containerul în acord cu regulamentele locale/regionale/naționale/internaționale.

(SE)

Varning

Kan skada fertiliteten eller det ofödda barnet. Orsakar allvarlig ögonirritation.

Använd skyddshandskar/skyddskläder/ögonskydd/ansiktsskydd. Vid bestående ögonirritation: Sök läkarhjälp. VID KONTAKT MED ÖGONEN: Skölj försiktigt med vatten i flera minuter. Ta ur eventuella kontaktlinser om det går lätt. Fortsätt att skölja. Innehållet / behållaren avfallshanteras enligt lokala / regionala / nationella / internationella föreskrifter.

(SI)

Pozor

Lahko škoduje plodnosti ali nerojenemu otroku. Povzroča hudo draženje oči.

Nositi zaščitne rokavice/zaščitno obleko/zaščitno za oči/ zaščitno za obraz. Če draženje oči ne preneha: poiščite zdravniško pomoč/oskrbo. PRI STIKU Z OČMI: previdno izpirajte z vodo nekaj minut. Odstranite kontaktno lečo, če jih imate in če to lahko storite brez težav. Nadaljujte z izpiranjem. Vsebino/vsebnik odstranite v skladu z lokalnimi/regionalnimi/narodnimi/mednarodnimi predpisi.

(SK)

Pozor

Môže spôsobiť poškodenie plodnosti alebo nenarodeného dieťaťa. Spôsobuje vážne podráždenie očí.

Noste ochranné rukavice/ochranný odev/ochranné okuliare/ochranu tváre. Ak podráždenie očí pretrváva: vyhľadajte lekársku pomoc/starostlivosť. PO ZASIAHNUTÍ OČÍ: Niekoľko minút ich opatrne vyplachujte vodou. Ak používate kontaktné šošovky a ak je to možné, odstráňte ich. Pokračujte vo vyplachovaní. Zneškodnenie obsahu/obalu v súlade s miestnymi/oblastnými/národnými/medzinárodnými nariadeniami.

Bio-Rad

3, boulevard Raymond Poincaré
92430 Marnes-la-Coquette France
Tel. : +33 (0) 1 47 95 60 00
Fax.: +33 (0) 1 47 41 91 33
www.bio-rad.com



2015/03
881177