

Acute Phase Response
 Cancer
 Cardiovascular Disease
 Cytokines, Chemokines,
 Growth Factors
 Diabetes
 Toxicology
 Genotyping
 Immunoglobulin Isotyping
 Signal Transduction

Bio-Plex Pro™ Human Chemokine Panel

6Ckine / CCL21, BCA-1 / CXCL13, CTACK / CCL27, ENA-78 / CXCL5, Eotaxin / CCL11, Eotaxin-2 / CCL24, Eotaxin-3 / CCL26, Fractalkine / CX3CL1, GCP-2 / CXCL6, GM-CSF, Gro- α / CXCL1, Gro- β / CXCL2, I-309 / CCL1, IFN- γ , IL-1 β , IL-2, IL-4, IL-6, IL-8 / CXCL8, IL-10, IL-16, IP-10 / CXCL10, I-TAC / CXCL11, MCP-1 / CCL2, MCP-2 / CCL8, MCP-3 / CCL7, MCP-4 / CCL13, MDC / CCL22, MIF, MIG / CXCL9, MIP-1 α / CCL3, MIP-1 δ / CCL15, MIP-3 α / CCL20, MIP-3 β / CCL19, MIPF-1 / CCL23, SCYB16 / CXCL16, SDF-1 α + β / CXCL12, TARC / CCL17, TECK / CCL25, TNF- α

MAGNETIC SEPARATION ENABLED

- All-in-one 40-plex kit
- Custom configurations
- Single level quality control
- Magnetic workflow



High-Performance Multiplex Immunoassays for Research

The Bio-Plex Pro Human Chemokine Panel features 40 magnetic bead-based immunoassays to measure chemokines and chemotaxis-relevant analytes. The assays have undergone rigorous evaluation to provide reliable performance for your research needs and are available in flexible, customized formats. Chemokines are typically associated with the following areas of research:

- Cancer/angiogenesis
- Cardiovascular disease
- Autoimmune disorders
- Allergy/asthma
- Diabetes, type 2
- Atherosclerosis
- COPD (chronic obstructive pulmonary disease)
- Dermatitis/psoriasis
- Alzheimer's disease

Assay Features

- Magnetic beads for simplified plate processing
- Single level quality control with kit lot-specific ranges
- Assay quick guide to get you started right away
- Compatible with Bio-Plex® 200, Bio-Plex 3D, and Bio-Plex® MAGPIX™ Systems

Rigorous Assay Validation

All Bio-Plex Pro Assays undergo rigorous evaluation that includes the following parameters:

- Specificity (cross-reactivity)
- Accuracy (recovery) in key sample matrices
- Inter- and intra-assay precision
- Sensitivity (limit of detection, LOD)
- Assay working range (LLOQ/ULOQ)
- Linearity of dilution
- Parallelism and matrix effect
- Performance characteristics in real samples

Assay Performance Definitions

The following parameters are indicative of assay performance, as shown in Table 1.

- **Assay working range** — the range of concentrations within which the assay is precise and accurate. Boundaries of the assay working range are defined by the lower limit of quantification (LLOQ) and the upper limit of quantification (ULOQ)
- **Precision** — the coefficient of variation (%CV) at concentrations within the assay working range
- **Accuracy (recovery)** — percentage of the observed concentration relative to the expected concentration of a known amount of analyte within the assay working range
- **Sensitivity (limit of detection, LOD)** — the concentration of analyte for which the fluorescence intensity signal is 2 standard deviations above the background signal

BIO-RAD

Table 1. Representative performance characteristics.

Analyte	Assay Working Ranges, pg/ml		Assay Sensitivity	Assay Precision		Calibration to WHO/NIBSC	
	LLOQ	ULOQ	LOD	Intra-Assay %CV	Inter-Assay %CV	Factor*	Catalog #
6Ckine / CCL21	21.9	3,923	12.0	4	6	–	–
BCA-1 / CXCL13	0.7	1,200	0.1	2	2	–	–
CTACK / CCL27	1.2	5,000	0.3	3	5	–	–
ENA-78 / CXCL5	7.3	120,000	5.7	3	6	–	–
Eotaxin / CCL11	1.5	3,859	0.7	3	4	–	–
Eotaxin-2 / CCL24	6.2	4,073	3.2	3	4	–	–
Eotaxin-3 / CCL26	0.9	12,109	0.5	3	4	–	–
Fractalkine / CX3CL1	4.0	11,463	0.9	3	4	–	–
GCP-2 / CXCL6	0.8	11,135	0.6	3	7	–	–
GM-CSF	5.3	35,000	1.0	4	3	–	–
Gro- α / CXCL1	3.1	7,024	4.2	2	4	–	–
Gro- β / CXCL2	4.6	13,257	2.7	3	8	–	–
I-309 / CCL1	1.8	1,015	1.6	3	6	–	–
IFN- γ	2.3	20,236	0.4	3	5	0.7	87/586
IL-1 β	0.4	7,000	0.1	2	4	1.1	86/680
IL-2	0.8	13,000	0.1	3	6	–	–
IL-4	1.2	4,804	1.0	2	4	1.2	88/656
IL-6	0.7	12,000	0.1	2	5	1.0	89/548
IL-8 / CXCL8	0.5	7,640	0.04	3	4	1.7	89/520
IL-10	1.3	18,708	0.9	3	3	1.8	93/722
IL-16	2.1	34,000	0.8	3	4	–	–
IP-10 / CXCL10	1.6	7,714	1.1	2	7	–	–
I-TAC / CXCL11	0.1	2,298	0.05	3	8	–	–
MCP-1 / CCL2	0.3	4,812	0.1	3	3	2.5	92/794
MCP-2 / CCL8	0.3	4,056	0.04	3	5	–	–
MCP-3 / CCL7	1.9	20,133	1.3	6	6	–	–
MCP-4 / CCL13	0.2	3,368	0.1	3	4	–	–
MDC / CCL22	0.9	14,649	0.5	3	3	–	–
MIF	23.1	377,724	15.4	2	7	–	–
MIG / CXCL9	1.8	19,600	1.1	3	7	–	–
MIP-1 α / CCL3	0.4	1,543	0.3	4	6	1.9	92/518
MIP-1 δ / CCL15	1.7	9,100	0.2	3	5	–	–
MIP-3 α / CCL20	0.3	4,675	0.1	2	3	–	–
MIP-3 β / CCL19	3.0	48,494	1.1	2	4	–	–
MPIF-1 / CCL23	1.0	14,450	0.5	3	3	–	–
SCYB16 / CXCL16	0.5	2,867	0.1	3	4	–	–
SDF-1 α + β / CXCL12	8.3	115,730	10.3	4	8	–	–
TARC / CCL17	1.7	430	1.1	4	5	–	–
TECK / CCL25	20.6	114,493	4.9	4	5	–	–
TNF- α	0.9	13,879	0.2	3	8	1.8	88/786

NIBSC, National Institute for Biological Standards and Control; WHO, World Health Organization.

The LLOQ, ULOQ, LOD, and inter-assay precision %CV are mean data determined from three independent multiplex assays in a serum-based matrix. Intra-assay %CV is derived from one representative assay. LLOQ and ULOQ are defined as the boundary standard curve points in which the performance specifications of individual standard points were met for a 10% intra-assay CV and recovery range of 70–130%. Data were generated using the magnetic workflow with the Bio-Plex Pro Wash Station.

* Factor x NIBSC value (pg/ml) = Bio-Plex value (pg/ml). Factors are based on average percentage recovery of the Bio-Plex Standards (in the linear range) relative to NIBSC standards run on the same assay plate.

Accuracy of Bio-Plex Pro Human Chemokine Panel

Linearity of dilution displays the ability of an assay to generate measured values from complex samples by comparing the accuracy over a range of sample dilutions. Bio-Plex Pro Assays are designed and validated with high linearity to return accurate results from complex matrices (Figure 1).

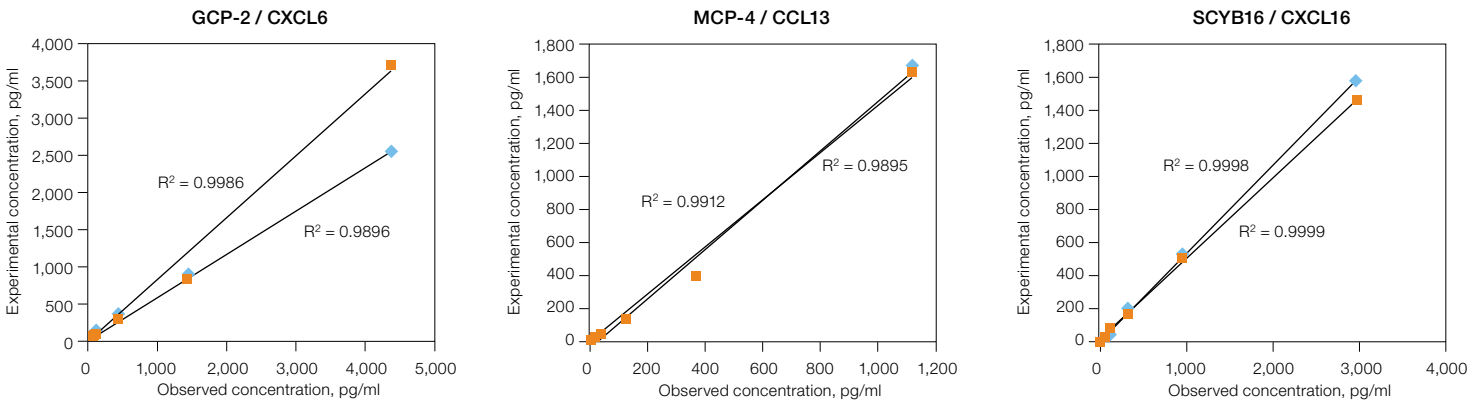


Fig. 1. Linearity of dilution determines the suitability of a standard curve for reflecting relative quantities of an analyte in a complex matrix. Linearity of dilution was assessed by spiking a known quantity of recombinant antigen into human serum and plasma matrices. Observed and expected analyte concentrations were plotted and the correlation coefficient (R^2) values reflect linearity in signal response. Plasma (◆); serum (■).

Bio-Plex Pro Assay Working Range

The assay working range should encompass the biological range of expression in order to be useful in research. Bio-Plex Pro Assays are developed and optimized to ensure real sample data fall within the quantifiable regions of the assay as demonstrated by comparing the standard curves of assay controls to biological samples (Figure 2).

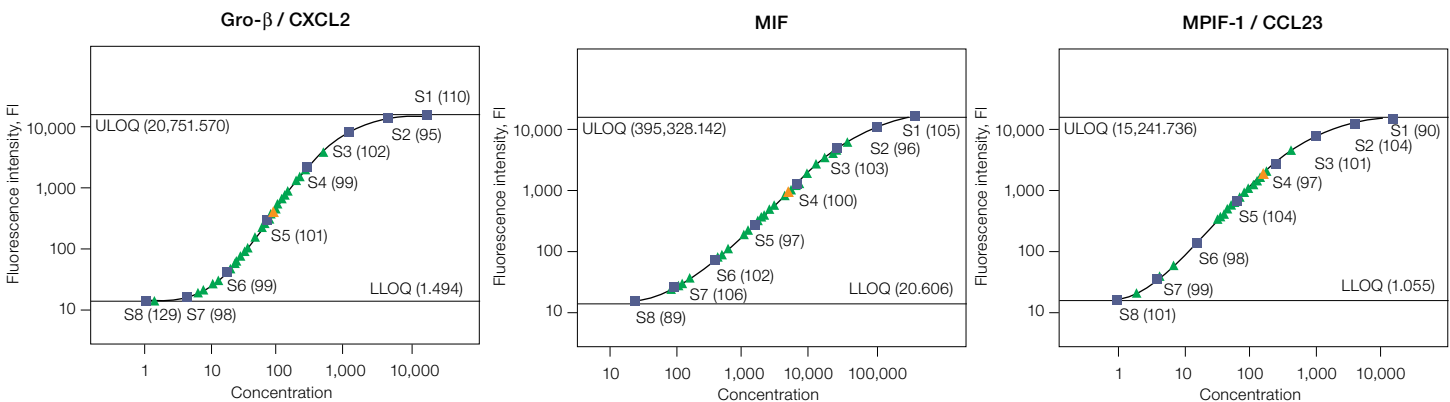


Fig. 2. Standard curves with assay controls and serum samples. Standard points were prepared by serially diluting a reconstituted standard fourfold to generate an eight-point standard curve. Standard points with percentage recovery (■); controls (▲); samples (▲). Data were generated in Bio-Plex Manager™ Software.

Detection of Analytes

Bio-Plex Pro Assays are tested with samples from multiple sources to ensure target analytes are detected within normal biological expression levels and levels associated with disease (Figure 3).

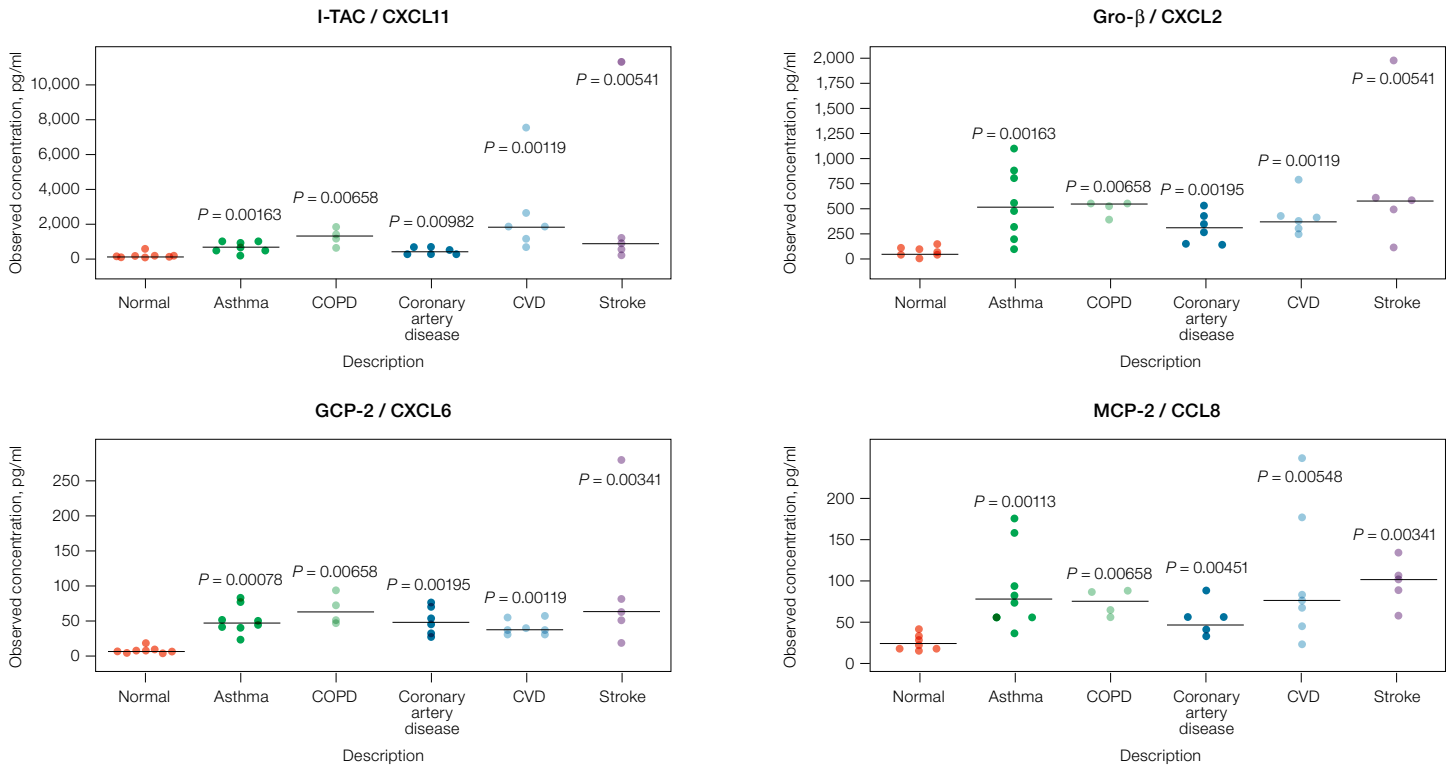


Fig. 3. Levels of biomarkers in sera of normal and various disease groups (asthma, COPD [chronic obstructive pulmonary disease], coronary artery disease, CVD [cardiovascular disease], and stroke). A Student *t* test was used to determine statistical significance between groups. Black lines denote mean values; *P* values are indicated above each population. Data analysis, graphing, and statistics were performed with Bio-Plex Data Pro™ Software.

Consistent Results across Different Multiplex Readers

Reliability of detection of analytes using Bio-Plex Pro Assays can be seen using a variety of Bio-Plex (or Luminex) platforms (Figure 4).

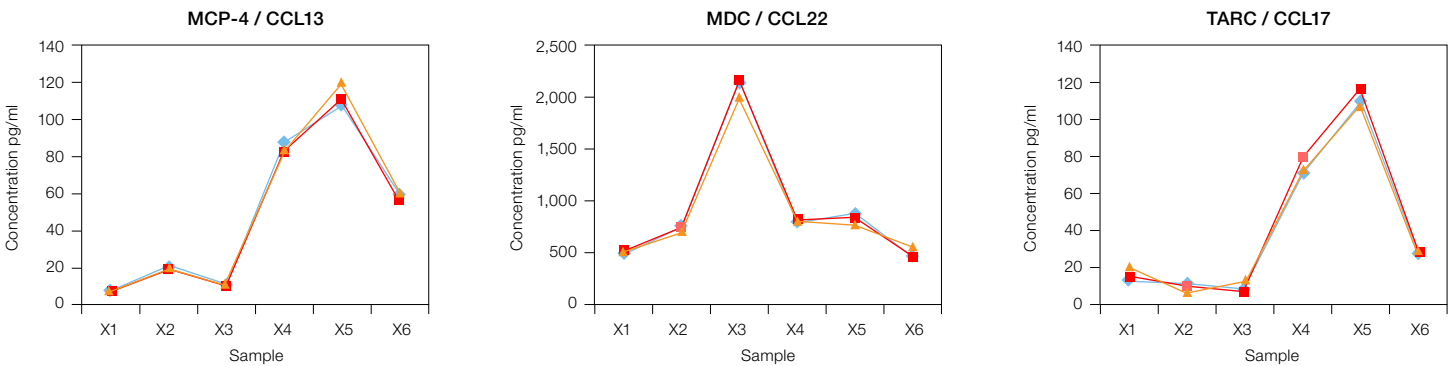


Fig. 4. Alignment of values from human serum samples across all Bio-Plex Readers. Bio-Plex MAGPIX (—▲—); Bio-Plex 200 (—■—); Bio-Plex 3D (—◆—). Data were generated in Bio-Plex Manager Software.

Flexible Ordering Options

Feature	Premixed Panel 40-Plex Panel	x-Plex™ Custom Assay (we mix)	Express Custom Assay (you mix)	Singleplex/Individual Components
All-in-one kit	•	•	•	
Single level quality controls	•			
Filter plate or flat bottom plate*	Flat bottom plate only	Filter or flat bottom plate	Filter or flat bottom plate	Filter or flat bottom plate
Choice of analytes		•	•	•
Number of analytes	40	11-plex or higher	10-plex or lower	Varies, 1–40
Detailed product data sheet	•	•	•	•
Assay quick guide**	•	•	•	•
Faster delivery	•		•	•
Premium performance	•	•		

* Filter plates are used for vacuum separation, flat bottom plates for magnetic separation.

** Go to bio-rad.com/web/Bio-PlexPanels for assay quick guides and manuals.

Ordering Information

Catalog #	Description	Catalog #	Description
Bio-Plex Pro Human Chemokines Panel All-in-One Kit		Reagent Kits	
171AK99MR2	Bio-Plex Pro Human Chemokine Panel , 1 x 96-well, includes coupled magnetic capture beads, premixed detection antibodies, standards, quality controls, detection antibody diluent HB, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions for the detection of 40 biomarkers noted below	171304090	Bio-Plex Pro Reagent Kit III with Filter Plate , 1 x 96-well, includes detection antibody diluent HB, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, filter plate, and sealing tape for vacuum separation methods
Bio-Plex Pro Human Chemokines Panel Singleplex Sets*		171304090M	Bio-Plex Pro Reagent Kit III with Flat Bottom Plate , 1 x 96-well, includes detection antibody diluent HB, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, flat bottom plate, and sealing tape for magnetic separation methods
171BK11MR2	6CKine / CCL21	Standards	
171BK12MR2	BCA-1 / CXCL13	171DK0001	Bio-Plex Pro Human Chemokine Standard , pkg of 1 vial, lyophilized mixture of 40 standard analytes
171BK13MR2	CTACK / CCL27	171DK0050	Bio-Plex Pro Human Chemokine Standard , pkg of 50 lot-matched vials, lyophilized mixture of 40 standard analytes
171BK14MR2	ENA-78 / CXCL5	Wash Stations and Accessories	
171BK15MR2	Eotaxin / CCL11	30034376	Bio-Plex Pro Wash Station , microplate wash station for magnetic bead-based assays, includes magnetic plate carrier, waste bottle, 2 liquid bottles
171BK16MR2	Eotaxin-2 / CCL24	171020100	Bio-Plex Handheld Magnetic Washer , includes magnetic washer and adjustment hex tools for use in manual wash steps for all Bio-Plex Magnetic Assays
171BK17MR2	Eotaxin-3 / CCL26	171025001	Bio-Plex Pro Flat Bottom Plates , pkg of 40, 96-well plates, for use with Bio-Plex Pro Wash Stations when using magnetic bead-based assays
171BK18MR2	Fractalkine / CX3CL1	Software	
171BK19MR2	GCP-2 / CXCL6	171001510	Bio-Plex Data Pro Software with Bio-Plex Manager Software , Bio-Plex Data Pro Software (5 seats), for multi-experiment analysis and advanced data visualization, and Bio-Plex Manager Software (5 seats), for instrument data evaluation and optimization. CDs and security HASP key included
171BK21MR2	GM-CSF	171001513	Bio-Plex Data Pro Software , (5 seats), for multi-experiment analysis and advanced data visualization
171BK22MR2	Gro- α / CXCL1	171STND01	Bio-Plex Manager Software , includes 1 user desktop license, for analysis of Bio-Plex data and generation of protocols, does not operate the instrument with Bio-Plex Pro Wash Stations when using magnetic bead-based assays
171BK23MR2	Gro- β / CXCL2	Visit bio-rad.com/web/Bio-PlexPanels for more information.	
171BK24MR2	I-309 / CCL1		
171BK25MR2	IFN- γ		
171BK26MR2	IL-1 β		
171BK27MR2	IL-2		
171BK28MR2	IL-4		
171BK29MR2	IL-6		
171BK31MR2	IL-8 / CXCL8		
171BK32MR2	IL-10		
171BK33MR2	IL-16		
171BK34MR2	IP-10 / CXCL10		
171BK35MR2	I-TAC / CXCL11		
171BK36MR2	MCP-1 / CCL2		
171BK37MR2	MCP-2 / CCL8		
171BK38MR2	MCP-3 / CCL7		
171BK39MR2	MCP-4 / CCL13		
171BK41MR2	MDC / CCL22		
171BK42MR2	MIF		
171BK43MR2	MIG / CXCL9		
171BK44MR2	MIP-1 α / CCL3		
171BK46MR2	MIP-1 δ / CCL15		
171BK47MR2	MIP-3 α / CCL20		
171BK48MR2	MIP-3 β / CCL19		
171BK49MR2	MPIF-1 / CCL23		
171BK51MR2	SCYB16 / CXCL16		
171BK52MR2	SDF-1 α + β / CXCL12		
171BK53MR2	TARC / CCL17		
171BK54MR2	TECK / CCL25		
171BK55MR2	TNF- α		

* Singleplex sets should not be mixed with others from different panels or groups.

HASP is a trademark of Aladdin Knowledge Systems, Ltd. Luminex is a trademark of Luminex Corporation.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666
New Zealand 64 9 415 2280 **Norway** 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

