

# Connectors for Using Bio-Rad and Cytiva Columns on NGC Chromatography Systems

## Quick Guide

### Columns and Recommended Connectors

Bio-Rad Laboratories, Inc. Columns							Cytiva Columns						
<b>EconoFit</b>	<b>Bio-Scale Mini</b>	<b>Foresight</b>	<b>UNO</b>	<b>Bio-Scale</b>	<b>ENrich</b>	<b>Econo Alpha</b>	<b>XK/HiLoad</b>	<b>Tricorn</b>	<b>HiScale</b>	<b>HiPrep</b>	<b>HiScreen</b>	<b>RESOURCE</b>	<b>HiTrap</b>

#### Recommended Bio-Rad Connectors

<b>A</b>	<b>B</b>	<b>C</b>	<b>D</b>	<b>E</b>	<b>F</b>
<b>10-32 PEEK 1/16"</b> (catalog #7800008)	<b>10-32 PEEK Union</b> (#7885007)	<b>Quick Connect Luer M to 10-32 F</b> (#7885010)	<b>Slip Luer F to 10-32 F</b> (#7885010)	<b>10-32 Narrow Fingertight 1/16"</b>	<b>1/4-28 Nut 1/16" and Ferrule 1/16"</b> (#7885025, 7500556)
Provided with NGC Systems and Column Switching Valves*	Included in NGC Fittings Kit (#7884017)	Included in NGC Fittings Kit (#7884017)	Included in NGC Fittings Kit (#7884017)	Provided with EconoFit Columns	Included in NGC Fittings Kit (#7884017)

\* Tefzel tubing with 1/4-28 fittings and ferrules provided with Bio-Scale Columns.

\*\* Econo Alpha Columns include a 1/4-28 M to 10-32 F adaptor.

**Bio-Rad and Cytiva Lab-Scale Chromatography Column Cross-Reference Guide**

Technique	Bio-Rad Columns	Comparable Cytiva Columns	Available Bio-Rad Column Formats
Affinity	EconoFit Nuvia IMAC, Ni-Charged	HiTrap FF, HiTrap HP	1 x 1 ml (#12009288), 5 x 1 ml (#12009285), 1 x 5 ml (#12009286), 5 x 5 ml (#12009287)
	EconoFit Nuvia IMAC, Uncharged	HiTrap IMAC FF, HiTrap IMAC HP	1 x 1 ml (#12009289)
	EconoFit Profinity IMAC, Ni-Charged	Capto Chelating*	5 x 1 ml (#12009298), 1 x 5 ml (#12009299), 5 x 5 ml (#12009300)
	EconoFit UNOsphere SUPrA	HiTrap rProtein A FF, HiTrap Protein A HP, HiTrap MabSelect	1 x 1 ml (#12009321), 5 x 1 ml (#12009322), 1 x 5 ml (#12009323)
	EconoFit Profinity GST**	GSTrap FF, GSTrap HP	5 x 1 ml (#12009295), 1 x 5 ml (#12009296), 5 x 5 ml (#12009297)
	EconoFit Affi-Gel Blue	HiTrap Blue HP	1 x 5 ml (#12009234), 5 x 5 ml (#12009235)
Ion Exchange	ENrich Q	Mono Q GL	1 x 1 ml (#7800001), 1 x 8 ml (#7800003)
	ENrich S	Mono S GL	1 x 1 ml (#7800021), 1 x 8 ml (#7800023)
	EconoFit UNOsphere Q	HiTrap Q FF	1 x 1 ml (#12009307), 5 x 1 ml (#12009301), 1 x 5 ml (#12009302), 5 x 5 ml (#12009303)
	EconoFit Macro-Prep High Q	HiTrap Q FF, HiTrap Q HP	1 x 1 ml (#12009275), 5 x 1 ml (#12009267), 1 x 5 ml (#12009268), 5 x 5 ml (#12009269)
	EconoFit UNOsphere S	HiTrap SP FF	1 x 1 ml (#12009308), 5 x 1 ml (#12009304), 1 x 5 ml (#12009305), 5 x 5 ml (#12009306)
	EconoFit Macro-Prep High S	HiTrap SP FF, HiTrap SP HP	1 x 1 ml (#12009276), 5 x 1 ml (#12009270), 1 x 5 ml (#12009271), 5 x 5 ml (#12009272)
	EconoFit Macro-Prep DEAE	HiTrap DEAE Sepharose FF	1 x 1 ml (#12009274), 5 x 1 ml (#12009264), 1 x 5 ml (#12009265), 5 x 5 ml (#12009266)
	EconoFit Macro-Prep CM	HiTrap CM FF	1 x 1 ml (#12009273)
	EconoFit Nuvia Q	HiTrap Q XL, HiTrap Capto Q	1 x 1 ml (#12009290)
	EconoFit Nuvia S	HiTrap SP XL, HiTrap Capto S	1 x 1 ml (#12009291)
	EconoFit Nuvia HP-Q	HiTrap Capto Q ImpRes	1 x 1 ml (#12009282)
	EconoFit Nuvia HR-S	HiTrap Capto SP ImpRes	1 x 1 ml (#12009284)
Mixed Mode	EconoFit CHT Type I***	HiTrap Capto adhere, HiTrap Capto adhere ImpRes	1 x 5 ml (#12009253), 5 x 5 ml (#12009254), 1 x 1 ml (#12009255)
	EconoFit CHT XT***	HiTrap Capto adhere, HiTrap Capto adhere ImpRes	1 x 1 ml (#12009261)
Hydrophobic Interaction	EconoFit Macro-Prep t-Butyl	HiTrap Butyl FF, HiTrap Butyl HP	1 x 1 ml (#12009278)
Size Exclusion/ Desalting	ENrich SEC 650	Superdex 200 GL, Superdex 200 Increase GL	1 x 24 ml (#7801650)
	ENrich SEC 70	Superdex 75 GL, Superdex 75 Increase GL	1 x 24 ml (#7801070)
	EconoFit Bio-Gel P-6 Desalting	HiTrap Desalting	1 x 5 ml (#12009238), 5 x 5 ml (#12009239)
	Bio-Scale Mini Bio-Gel P-6 Desalting	HiPrep 26/10 Desalting	1 x 10 ml (#7325304), 1 x 50 ml (#7325312), 5 x 50 ml (#7325314)
	Econo-Pac 10DG Desalting	PD-10 Desalting	30 x 10 ml (#7322010)

\* Not available in column format.

\*\* Consider also EconoFit Profinity eXact Columns, which enable tag-free protein purifications.

\*\*\* CHT Ceramic Hydroxyapatite Media and Capto adhere/adhere ImpRes Media are chemically distinct but exhibit superior performance for monoclonal antibody purification workflows (see [bulletin 6749](#)). CHT Media offer unique selectivities and often separate biomolecules that other chromatography supports are unable to differentiate.Visit [bio-rad.com/Columns](http://bio-rad.com/Columns) to view our full offering of Chromatography Columns.Visit [bio-rad.com/EconoFit](http://bio-rad.com/EconoFit) to view our full selection of EconoFit Prepacked Columns.Visit [bio-rad.com/NGC](http://bio-rad.com/NGC) for more information on our NGC Chromatography Systems.

BIO-RAD, AFFI-GEL, BIO-GEL, BIO-SCALE, CHT, ECONO-PAC, MACRO-PREP, and UNO are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

**BIO-RAD****Bio-Rad  
Laboratories, Inc.**Life Science  
Group

**Website** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050  
**Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670  
**The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23  
**Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23  
**Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

