

HIS-Select Quick Reference Guide



	Nickel Affinity Gel	High Flow (HF) Nickel Affinity Gel	Cobalt Affinity Gel	Nickel Cartridges	Nickel Spin Columns	EZview™ Red Affinity Gel	Nickel Magnetic Beads	High Sensitivity (HS) Nickel Coated Strip Plates	High Capacity (HC) Nickel Coated Plates 96-well	iLAP™ High Capacity (HC) Nickel Coated Plates 96-well	Filter Plates 96-well
Product Number	P 6611	H 0537	H 8162	H 8286	H 7787	E 3528	H 1786	S 5688	S 5563, H 1661	H 9412	Coming Fall 2003
Package Sizes	5 ml, 25 ml, 100 ml, 500 ml	10 ml, 25 ml, 100 ml, 500 ml	5 ml, 25 ml, 100 ml	5 ea	10 ea, 50 ea	1 ml, 5 x 1 ml	2 x 1 ml, 5 x 1 ml	1 ea, 5 ea	1 ea, 5 x 1 ea	1 ea, 5 x 1 ea	1 ea, 5 ea
Application	<ul style="list-style-type: none"> • Gravity Flow Column • Small-Med. Scale 	<ul style="list-style-type: none"> • FPLC™ • Production Scale 	<ul style="list-style-type: none"> • Gravity Flow Column • Small-Med. Scale 	<ul style="list-style-type: none"> • FPLC • Syringe Purification • Small Scale 	<ul style="list-style-type: none"> • Mini Prep • Process by Centrifugation or Vacuum 	<ul style="list-style-type: none"> • Mini Prep • Pull-Downs 	<ul style="list-style-type: none"> • Rapid Mini Prep Magnetic Pull-Downs • High Throughput Screening • Compatible with Robotics when used in 96-well plates 	<ul style="list-style-type: none"> • High Throughput Screening (ELISA) • Compatible with Robotics 	<ul style="list-style-type: none"> • High Throughput, High Capacity Screening • Multi-Sample Mini Prep • In-Well Protein Assays • Compatible with Robotics 	<ul style="list-style-type: none"> • Direct Cell Lysis and Purification of Target Protein in Well • High Throughput, High Capacity Screening • Multi-Sample Mini Prep • In-Well Protein Assays • Compatible with Robotics 	<ul style="list-style-type: none"> • Small Scale • High Throughput Screening • Process by Centrifugation or Vacuum • Compatible with Robotics
Scale per Sample	100 µg – 10 g	100 µg – 100 g	100 µg – 10 g	15 – 20 mg	<ul style="list-style-type: none"> • 150 µg • 600 µl/load 	10 – 500 µg per pull-down	10 µg – 15 mg per pull-down	<ul style="list-style-type: none"> • 1 ng – 1 µg/well • 96-well: 200 µl/well 	<ul style="list-style-type: none"> • 1 – 4 µg/well • 96-well: 200 µl/well • 384-well: 60 µl/well 	<ul style="list-style-type: none"> • 1 – 4 µg/well • 96-well: 200 µl/well 	<ul style="list-style-type: none"> • 1 mg/well • 1 ml/load
Binding Capacity	15 mg/ml	15 mg/ml	15 mg/ml	15 mg/cartridge	150 µg/column	15 mg/ml	15 mg/ml	Sensitivity: ≤1 ng/well	96-well: ≥4 µg/well 384-well: ≥ 2µg/well	≥4 µg/well	1 mg/well
Matrix	6% Beaded Agarose	6% Beaded Agarose, Highly Cross-linked	6% Beaded Agarose	6% Beaded Agarose	Silica Particles (Porous, Spherical)	6% Beaded Agarose	Paramagnetic Iron Impregnated 6% Beaded Agarose	Clear polystyrene plate; Proprietary Coating	Clear Polystyrene Plate; Proprietary Coating	Clear polystyrene plate; Proprietary Coatings for lysis and purification	Silica Particles (Porous, Spherical)
Bead Size	45 – 165 µm	45 – 165 µm	45 – 165 µm	45 – 165 µm	Approx. 20 µm	45 – 165 µm	20 – 75 µm	N/A	N/A	N/A	Approx. 20 µm
Exclusion Limit (MW) or Pore Size	4 x 10 ⁶ Da	4 x 10 ⁶ Da	4 x 10 ⁶ Da	4 x 10 ⁶ Da	1000 Angstrom (approx. 10 x 10 ⁶ Da)	4 x 10 ⁶ Da	4 x 10 ⁶ Da	N/A	At least 300 kDa	At least 300 kDa	1000 Angstrom (approx. 10 x 10 ⁶ Da)
Max. Linear Flow Rate (Max. pressure)	150 cm/hr (5 psi)	3000 cm/hr (200 psi)	150 cm/hr (5 psi)	150 cm/hr (5 psi)	N/A	N/A	N/A	N/A	N/A	N/A	N/A
Recommended Flow Rate for 1 x 2 cm Column	1 ml/min	5 ml/min	1 ml/min	1 ml/min	N/A	N/A	N/A	N/A	N/A	N/A	N/A
Recommended Binding Time/Speed	N/A	N/A	N/A	N/A	Equivalent to approx. 80 x g Centrifugation	1 hr @ 4 °C	0.5 hr @ 25 °C	1 hr @ 25 °C	96-well: 4 hr @ 25 °C 384-well: 2 hr @ 25 °C	1-4 hr @ 25 °C	Equivalent to approx. 270 x g Centrifugation
Optimal pH Stability	3-10	3-10	3-10	3-10	3-10	3-10	3-10	3-10	3-10	3-10	3-10
Physical Form	50% suspension in 30% ethanol	50% suspension in 30% ethanol	50% suspension in 30% ethanol	1.25 ml pre-packed cartridge	Pre-packed dry matrix in spin column	50% suspension in 30% ethanol	50% suspension in 30% ethanol	Clear 96-well-flat bottom, polystyrene strip plate	Clear 96- and 384-well flat bottom, polystyrene plates	Clear 96-well flat bottom, polystyrene plates	Pre-packed dry matrix in 96-well filter plate
Antimicrobial Agent	30% ethanol	30% ethanol	30% ethanol	30% ethanol	N/A	30% ethanol	30% ethanol	N/A	chlorhexidine	N/A	N/A
Storage	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C	2-8 °C
Recommended Imidazole Conc. for Load/Wash	0-10 mM	0-10 mM	0-10 mM	0-10 mM	0-5 mM	0-10 mM	0-10 mM	0-5 mM	0-1 mM	0-1 mM	0-5 mM
Recommended Imidazole Conc. for Elution	250 mM	250 mM	250 mM	250 mM	250 mM	250 mM	250 mM	N/A	200 mM	200 mM	250 mM

