

# 蛋白质纯化工具

从表达到检测, 帮助您获得更佳结果

应用	靶点	配体/化学特性	Thermo Scientific® 和 Invitrogen® 产品	磁珠/树脂类型	结合能力 (详情请参见网站)	最高流速 (cm/hr)	包装 (参见图1)	推荐规模 (参见表1)				
离子交换纯化	蛋白质电荷	强阴离子交换	thermofisher.com/sax	POROS® HQ 树脂 耐盐性 = 150 mM	60-70 mg/mL (动态载量)	5,000	B	B	P	PRO		
			thermofisher.com/scx	POROS® XQ 树脂	>140 mg/mL (动态载量)			B	P	PRO		
			thermofisher.com/ionexchange	POROS® XS 树脂	>100 mg/mL (动态载量)			B	P	PRO		
融合蛋白纯化	His 标签蛋白	Ni-NTA (更高的结合能力)	thermofisher.com/ninta	HisPur® Ni-NTA 磁珠	≥40 µg/mg磁珠	不适用	HTB	B				
			thermofisher.com/ninta	Pierce® Ni-NTA 磁性琼脂糖	>70 mg/mL	不适用		B				
			thermofisher.com/ninta	HisPur® Ni-NTA 树脂	~60 mg/mL	700		B				
			thermofisher.com/ninta	HisPur® Ni-NTA Superflow 琼脂糖	>60 mg/mL (动态载量)	1,200		B	P			
	钴 (更高的纯度)	HisPur® 钴树脂	thermofisher.com/cobalt	HisPur® 钴树脂	≥15 mg/mL	700	B					
			thermofisher.com/cobalt	HisPur® 钴Superflow琼脂糖	>30 mg/mL (动态载量)	1,200	B	P				
	GST 融合蛋白	谷胱甘肽	thermofisher.com/gstpuration	Pierce® 谷胱甘肽磁性琼脂糖	>12 mg/mL	不适用	HTB	B				
				Pierce® 谷胱甘肽琼脂糖	~40 mg/mL	700		B				
				Pierce® 谷胱甘肽Superflow 琼脂糖	~30 mg/mL (动态载量)	1,200		B	P			
	DYKDDDDK(Flag) 标签蛋白	抗DYKDDDDK(Flag)抗体	thermofisher.com/flag	Pierce™ 抗DYKDDDDK磁性琼脂糖	≥3.2 mg/mL	NA	HTB	B				
				Pierce™ 抗DYKDDDDK聚丙烯酰胺亲和树脂	≥3 mg/mL	150		B	P			
	HA 标签蛋白	抗HA抗体	thermofisher.com/ha	Pierce® 抗HA磁珠	≥100 µg/mL bead suspension	NA	HTS	B				
Pierce® 抗HA琼脂糖				102-144 nmol/mL	700	B						
c-Myc 标签蛋白	抗c-Myc抗体	thermofisher.com/myc	Pierce® 抗c-Myc磁珠	≥100 µg/mL bead suspension	NA	HTS	B					
			Pierce® 抗c-Myc琼脂糖	60-150 nmol/mL	700		B	P				
抗体纯化	IgG	抗体	阴性筛选	Melon® 凝胶	大多数抗体可获得>90%的产量和>80%的纯度	700	B					
			蛋白 A	thermofisher.com/protein-a	Dynabeads® 蛋白A磁珠	25-30 µg/mg磁珠	不适用	HTS	B			
				thermofisher.com/protein-a	Pierce® 蛋白A Plus琼脂糖	34 mg/mL	700		B			
				thermofisher.com/protein-a	POROS® MabCapture® A Select 树脂	≥37 mg/mL (动态载量)	5,000		B	P	PRO	
			蛋白 G	thermofisher.com/protein-g	Dynabeads® 蛋白G磁珠	25-30 µg/mg磁珠	NA	HTS	B			
				thermofisher.com/protein-g	Pierce® 蛋白G琼脂糖	11-15 mg/mL	700		B			
				thermofisher.com/protein-g	POROS® MabCapture® G Select 树脂	≥13 mg/mL (动态载量)	5,000		B	P	PRO	
			蛋白 A/G	thermofisher.com/protein-ag	Pierce® 蛋白A/G磁珠	≥55-85 µg/mg磁珠	不适用	HTS	B			
				thermofisher.com/protein-ag	Pierce® 蛋白A/G磁性琼脂糖	>40 mg/mL			HTB			
				thermofisher.com/protein-ag	Pierce® 蛋白A/G琼脂糖	≥7 mg/mL			700	B		
				thermofisher.com/protein-ag	Pierce® 蛋白A/G Plus琼脂糖	≥50 mg/mL				B		
				thermofisher.com/protein-ag	POROS® MabCapture® A/G Select 树脂	≥22 mg/mL (动态载量)				5,000	B	P
			蛋白 L	thermofisher.com/protein-l	Pierce® Protein L 磁珠	≥110 µg/mg磁珠	不适用	HTS	B			
				thermofisher.com/protein-l	Pierce® 蛋白L Plus琼脂糖	5-10 mg/mL			B			
				thermofisher.com/protein-l	Pierce® 蛋白L琼脂糖	10-20 mg/mL			B			
嗜硫吸附剂	Pierce® 嗜硫吸附剂	thermofisher.com/sulf	Pierce® 嗜硫吸附剂	20 mg/mL	700	B						
		thermofisher.com/sulf	Pierce® 嗜硫吸附剂	20 mg/mL	700	B						
IgM	固相化甘露聚糖结合蛋白 (MBP)	thermofisher.com/mbp	Pierce® 甘露聚糖结合蛋白琼脂糖	1 mg/mL	700	B						
			Pierce® 甘露聚糖结合蛋白琼脂糖	1 mg/mL	700	B						
IgA	固相化葡聚糖	thermofisher.com/iga	Pierce® 葡聚糖琼脂糖	1-3 mg/mL	700	B						
			Pierce® 葡聚糖琼脂糖	1-3 mg/mL	700	B						
生物素亲和纯化	生物素化蛋白	生物素	亲和素	Pierce® 亲和素琼脂糖树脂	≥20 µg生物素/mL	700	B					
			亲和素	Pierce® 亲和素琼脂糖树脂	≥1.2 mg生物素化蛋白/mL	700	B					
			链霉亲和素	Pierce® 链霉亲和素磁珠	~55 µg生物素化蛋白/mg磁珠	不适用	HTS	B				
			链霉亲和素	Pierce® 高容量链霉亲和素琼脂糖	≥10 mg生物素化蛋白/mL	700		B				
链霉亲和素	NeutrAvidin® 中性亲和素蛋白	Pierce® NeutrAvidin® 中性亲和素琼脂糖	≥1-2 mg生物素化蛋白/mL	700	B							
蛋白质固定化	可与特定配基共价结合的蛋白	与蛋白氨基反应	Pierce® NHS 活化琼脂糖	>25 mg/mL	700	B						
			与蛋白巯基反应	AminoLink® 加强型偶联树脂	15-20 mg/mL	700	B					
			与蛋白巯基反应	SulfoLink® 偶联树脂	20 mg/mL	700	B					

表1. 根据纯化规模和应用选择树脂。

规模	HTS (高通量筛选)	HTB (高通量批量)	B (批量)	P (中试)	PRO (生产)
具体描述	小规模 自动化兼容性	实验室规模	实验室规模	根据要求扩大规模	生产规模
产量	毫克级	毫克级	毫克级	毫克级	毫克级
形式	磁珠处理系统	磁珠处理系统 多孔离心板 (琼脂糖)	重力流 离心柱 (琼脂糖)	中等流速FPLC	高速流FPLC
应用	高通量筛选 相互作用研究 突变分析	需要毫克级的高通量筛选	功能分析 结构分析	结构分析	大规模生产
推荐的树脂类型	磁珠 (1-2.8 µm)	磁性琼脂糖 (10-40 µm)	琼脂糖 (45-165 µm)	Superflow (45-165 µm)	UltraLink® resin (50-80 µm) POROS® 树脂 (50 µm)



图1. 包装形式概述。