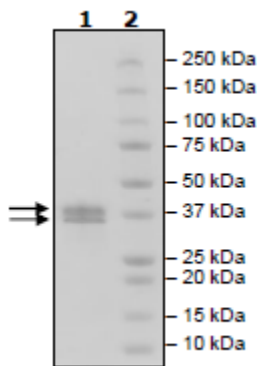


Product Information

Construct:	HMGB1 (2-215(end)-Avi-(biotin)-His)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%
Concentration:	0.45 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_002128
MW:	25 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

