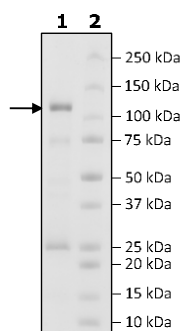


Product Information

Construct:	MDM2 (GST-Pre-2-498(end))
Concentration:	0.24 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT
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_002392
MW:	83 kDa
Purity:	52%
Assay Conditions:	Assay was performed according to MDM2 TR-FRET Assay Kit (BPS Bioscience #79773) with MDM2 titrated from 157.5 ng- 0 ng.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



MDM2 Activity

