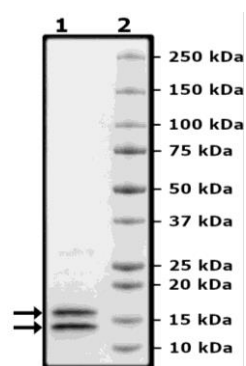


Product Information

Construct:	Caspase3 (1-277(end)-His)
Concentration:	0.90 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, and 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_004346
MW:	Small active subunit: 13 kDa; Large active subunit: 17 kDa
Purity:	≥90%
Specific Activity:	6784 pmol/min/g
Assay Conditions:	Assay was done according to Caspase-3 Homogeneous Assay Kit (BPS Bioscience #80700) with various amounts of Caspase-3.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Caspase3 Activity

