

Product Information

Description:	Recombinant human Dicer (Human Dicer, also known as Dicer 1, Ribonuclease III), full length, encompassing amino acids 2- 1922(end). This construct contains an N-terminal His-Tag (6xHis). This recombinant protein was affinity purified.
Species	Human
Construct:	Dicer (His-2-1922(end))
Concentration:	0.20 mg/ml
Expression System:	Sf9
Purity:	55%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, and 20% glycerol
MW:	221 kDa
Genbank Accession:	NM_030621
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

