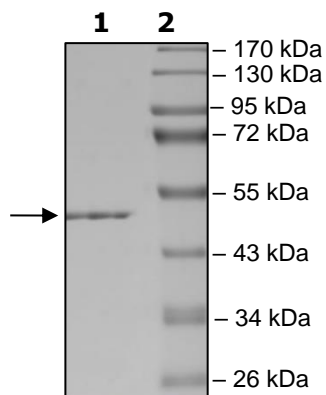


## Product Information

<b>Description:</b>	Recombinant <i>T. thermophilus</i> TtCsm6 (type III-A CRISPR-associated protein Csm6). This construct contains a C-terminal His-tag. TtCsm6 is a single-stranded RNA-specific endoribonuclease protein that contains an N-terminal CARF domain and C-terminal HEPN domain.
<b>Species:</b>	<i>Thermus thermophilus</i>
<b>Construct:</b>	Csm6 (Full Length-His) ( <i>T. thermophilus</i> )
<b>Concentration:</b>	0.20 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM sodium phosphate, pH 7.5, 300 mM NaCl, 1 mM DTT, 150 mM imidazole, and 10% glycerol
<b>MW:</b>	53 kDa
<b>Genbank Accession:</b>	WP_011229148
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Assay Conditions:</b>	The TtCsm6 RNase activity was determined using a fluorogenic substrate, with fluorescent signal measured over time.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



TtCsm6 RNase Activity

