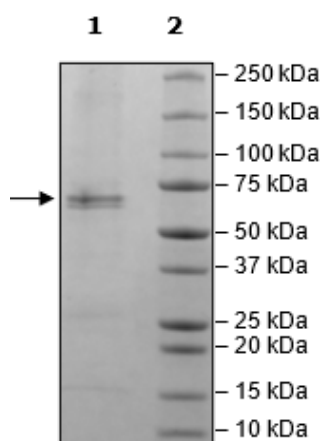


## Product Information

<b>Description:</b>	Recombinant Human PRMT1 (protein arginine N-methyltransferase 1), encompassing a.a. 2- 371 (end). This construct contains an N-terminal GST-tag followed by a thrombin cleavage site. The protein was affinity purified.
<b>Species:</b>	Human
<b>Construct:</b>	PRMT1 (GST-Th-2-371(end))
<b>Concentration:</b>	0.19 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	81%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT
<b>MW:</b>	69 kDa
<b>Genbank Accession:</b>	NM_001536
<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>	Assay was performed according to PRMT1 Chemiluminescent Assay Kit (BPS Bioscience #52004L) with various amounts of PRMT1.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PRMT1 Activity

