

## Product Information

<b>Description:</b>	Human Protein Arginine Methyltransferase 9 (PRMT9), full length encompassing amino acids 2-845 (end). This construct contains an N-terminal FLAG-tag and a C-terminal His-tag (6 x His). The recombinant protein was affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	PRMT9 (FLAG-2-845(end)-His)
<b>Concentration:</b>	2.16 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	≥90%
<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, 80 µg/ml FLAG peptide, and 2 mM TCEP
<b>MW:</b>	95 kDa
<b>Genbank Accession:</b>	NM_138364
<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.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

