

## Product Information

<b>Description:</b>	Recombinant human SET8 (PR/SET domain containing protein 8), encompassing amino acids 195-352. This construct contains an N-terminal GST-tag. This protein was affinity purified
<b>Background:</b>	SET8 (PR/SET domain containing protein 8), also known as KMT5A (lysine-specific methylates 5A), belongs to the family of proteins that contain an SET domain. It has the specificity of being the only protein that monomethylates H4K20me1, but also participates in monomethylating proteins like p53 and PCNA (proliferating cell nuclear antigen). It is involved in cell cycle regulation and DNA replication and repair. Abnormal expression of SET8 can lead to cancer, and this protein is found at high levels in several cancers, such as breast cancer. Inhibition of SET8 was found to increase cancer cell sensitivity to radiotherapy. The use of inhibitors alone or in combinatory therapy promises to bring benefits in cancer therapy.
<b>Species:</b>	Human
<b>Construct:</b>	Set8 (GST-195-352(end))
<b>Concentration:</b>	1.81 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<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, 20% glycerol, 3 mM DTT, and 3 mM Glutathione
<b>MW:</b>	44 kDa
<b>Genbank Accession:</b>	NM_020382
<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>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-PAGE Coomassie Staining

