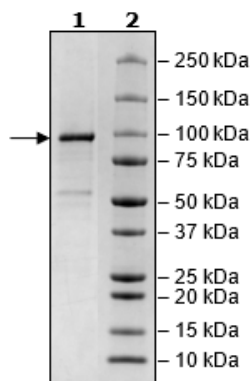


## Product Information

<b>Construct:</b>	PDE3B (GST-592-1112(end))
<b>Concentration:</b>	0.45 mg/ml
<b>Species:</b>	Human
<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>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_000922
<b>MW:</b>	87 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	1270 pmol/min/μg
<b>Assay Conditions:</b>	Assay was performed with 10 mM Tris-HCl, pH 7.4, 10 mM MgCl <sub>2</sub> , 0.05% Tween-20, 200 μM cGMP, 2.5 kU 5'-nucleotidase, 0.1 mg/ml BSA, and material serially diluted. The reaction was incubated at 37°C for 20 min. Following the reaction, malachite green reagent was added and then read 10 min after with a machine capable of reading absorbance at 620 nm.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



PDE3B Activity

