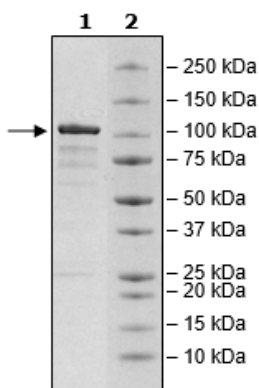


Product Information

Construct:	PDE4B1 (GST-1-736(end))
Concentration:	0.71 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, and 20% glycerol
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001037341
MW:	109 kDa
Purity:	81%
Specific Activity:	8532 pmol/min/μg
Assay Conditions:	Assay was performed with PDE Assay Buffer (BPS Bioscience #60393), 0.1 mg/ml BSA, 500 μM cAMP, 0.05 Ku/μl 5' Nucleotidase (Enzo-BML-K1307), and PDE4B1, GST-Tag (BPS Bioscience #60041) serial diluted at various concentrations. 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 microplate reader capable of reading absorbance at 620 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



PDE4B1 Activity

