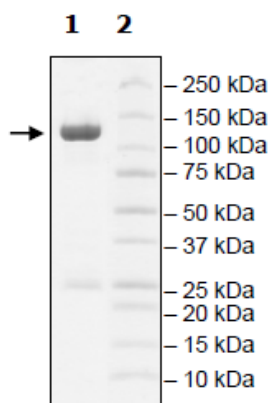


Product Information

Construct: PDE11A4 (GST-1-934(end))
Concentration: 0.93 mg/ml
Species: Human
Formulated In: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 3 mM DTT, 1.6 mM glutathione
Expression System: Sf9
Format: Aqueous buffer solution
Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage: -80°C
Genbank Accession: BC112393
MW: 131 kDa
Purity: ≥90%
Specific Activity: 1560 pmol/min/μg
Assay Conditions: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.1 mg/ml BSA, 0.05% Tween-20, 200 μM cAMP, and serial dilutions of PDE11A4 at 37°C for 20 min. Quantified by 5' nucleotidase cleaving the 5'-AMP product and releasing the phosphate group which is detected by Malachite Green Reagent.
Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

