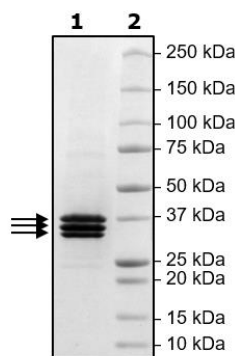


Product Information

Construct:	CDK7 (His-2-346(end)) / CyclinH1 (1-323(end)-His) / MNAT1 (His-2-309(end))
Concentration:	1.09 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 117 mM imidazole
Expression System:	Coexpressed in Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	CDK7: NM_001799; CyclinH1: NM_001239; MNAT1: NM_002431
MW:	CDK7: 40 kDa; CyclinH1: 39 kDa; MNAT1: 37 kDa
Purity:	≥90%
Specific Activity:	21 pmol/min/μg
Assay Conditions:	Assay was done according to the CDK7 Assay Kit (BPS Bioscience #79603) with CDK7/CyclinH1/MNAT1 titrated from 40 ng/rxn to 0.625 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

