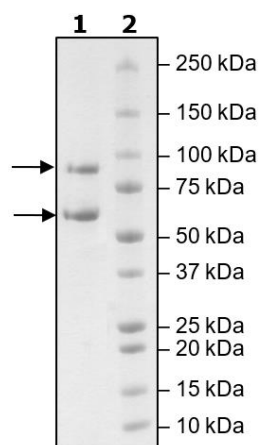


Product Information

| | |
|------------------------------|---|
| Description: | Complex of human CDK1 (cyclin-dependent kinase 1) and recombinant human cyclin B1 (CCNB1). The proteins were co-expressed in Sf9 insect cells. Recombinant human CDK1, full length, encompassing amino acids 2-297(end). This construct contains an N-terminal GST-tag followed by a PreScission Protease Cleavage Site. Recombinant human cyclin B1, full length, encompassing amino acids 2-433(end). This construct also contains an N-terminal GST-tag followed by a PreScission Protease Cleavage Site. The recombinant proteins were affinity purified. |
| Species | Human |
| Construct: | CDK1 (GST-297(end)) / CyclinB1 (GST-2-433(end)) |
| Concentration: | 0.50 mg/ml |
| Expression System: | Sf9 |
| Purity: | ≥90% |
| Format: | Aqueous buffer solution. |
| Formulated In: | 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT |
| MW: | CDK1: 60 kDa; CyclinB1: 75 kDa |
| Genbank Accession: | CDK1: NM_001786; CyclinB1: NM_031966 |
| Stability: | At least 6 months at -80°C. |
| Storage: | -80°C |
| Instructions for Use: | 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. |
| Specific Activity: | ≥200 pmol/min/μg |
| Assay Conditions: | Assay was done according to CDK1 Assay Kit (BPS Bioscience #79597) with various amounts of CDK1. |
| Applications: | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CDK1/CyclinB1 Activity

