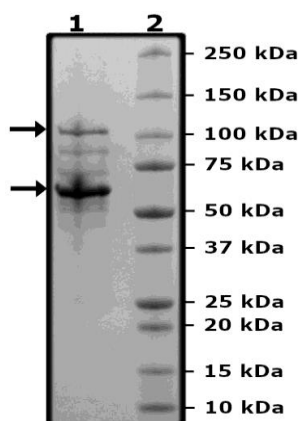


Product Information

| | |
|------------------------------|---|
| Description: | Complex of recombinant human cyclin dependent kinase 9 (CDK9), also known as cell division protein kinase 9, and recombinant human cyclin T1 (CCNT1, CycT1). The CDK9 construct is full length, encompassing amino acids 2-372(end), with N-terminal GST-tag. The cyclin T1 construct is full length, encompassing amino acids 2-726(end), with N-terminal GST-tag. The recombinant proteins were co-expressed and affinity purified. |
| Species | Human |
| Construct: | CDK9 (GST-2-372(end)) / CyclinT1 (GST-2-726(end)) |
| Concentration: | 0.11 mg/ml |
| Expression System: | Sf9 |
| Purity: | 77% |
| Format: | Aqueous buffer solution. |
| Formulated In: | 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.02% Tween-20, and 20% glycerol |
| MW: | CDK9: 69 kDa; CyclinT1: 107 kDa |
| Genbank Accession: | CDK9: NM_001261; CyclinT1: NM_001240 |
| 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: | 26 pmol/min/μg |
| Assay Conditions: | Assay was done according to CDK9 Assay Kit (BPS Bioscience #79628) with various amounts of CDK9. |
| Applications: | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CDK9/CyclinT1 Activity

