

Data Sheet

CDK7/CyclinH1/MNAT1, His-tags

Human, Recombinant, N-terminal His-Tags
Catalog #: 40098
Lot #: 140915 **Conc.:** 0.744 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04%
 Tween20, 20% glycerol, and 226mM
 imidazole

Stability: 12 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

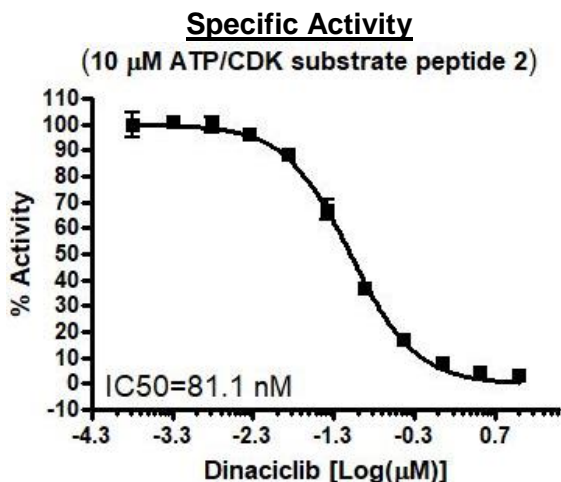
1. Fisher, R. P.; *Cell* **78**: 713-724, 1994.
2. Larochelle, S. *Molec. Cell* **25**: 839-850, 2007.

Description: Complex of recombinant human CDK7 (a.a. 2-346 [end]), Cyclin H1 (a.a. 1-323 [end]), and MNAT1 (a.a. 2-309 [end]) with N-terminal His-tags (CDK7 and MNAT1) and C-terminal His-tag (CyclinH1), GenBank Accession Nos. NM_001799, NM_001239, and NM_002431 respectively, co-expressed in Sf9 insect cells via a baculovirus expression system.

Assay Conditions: Inhibition of CDK7/Cyclin H/MAT1 enzyme by Dinaciclib, measured using the CDK7 assay kit (Cat. #79603).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 CDK7/cycH/MNAT
 (7 μg loaded)
Lane 2:
 Protein Marker

Purity: $\geq 60\%$

MW:
 CDK7: 40 kDa
 CyclinH: 38.5 kDa
 MNAT1: 37 kDa

