

CTK/MATK Kinase Assay Service

Activity-based CTK/MATK kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

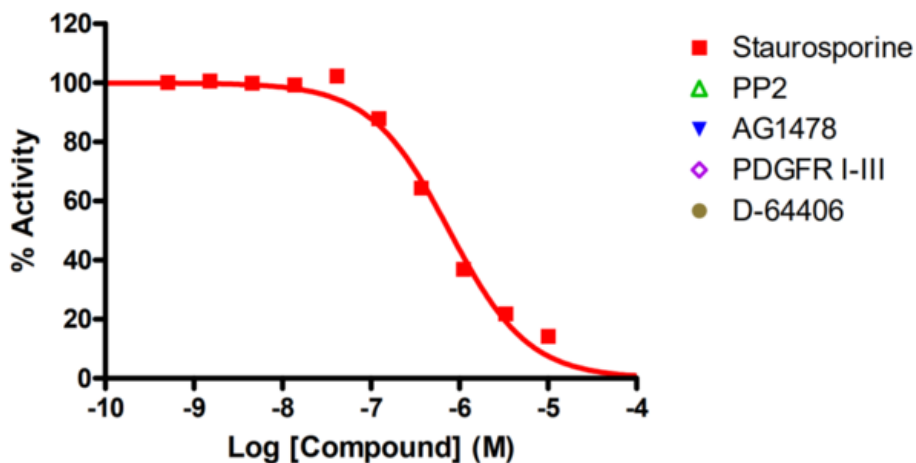
Target Overview

| | |
|------------|--|
| CAT# | CTK |
| Enzyme | Human CTK |
| HGNC | MATK |
| Synonym | CHK, CTK, DKFZp434N1212, HHYLTk, HYL, Lsk, MGC1708, MGC2101 |
| Substrates | Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml ATP 10 μM |

Assay Properties

| | |
|--------------------------|--|
| Reaction | Substrate + [³² P]-ATP = ³² P-Substrate + ADP |
| Readout | Enzymatic activity |
| Reference Compound IC50s | Compounds IC50 (nM) Staurosporine 770 |
| Scientific Information | P42679 (http://www.uniprot.org/uniprot/P42679) |
| Screening Location | Malvern, USA |

Reference compound IC50 for CTK/MATK/HYL



Let's Discover Together.

© Aug 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/ctk_matk_kin_malvern

