

BTK (E41K) Kinase Assay Service

Activity-based BTK (E41K) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

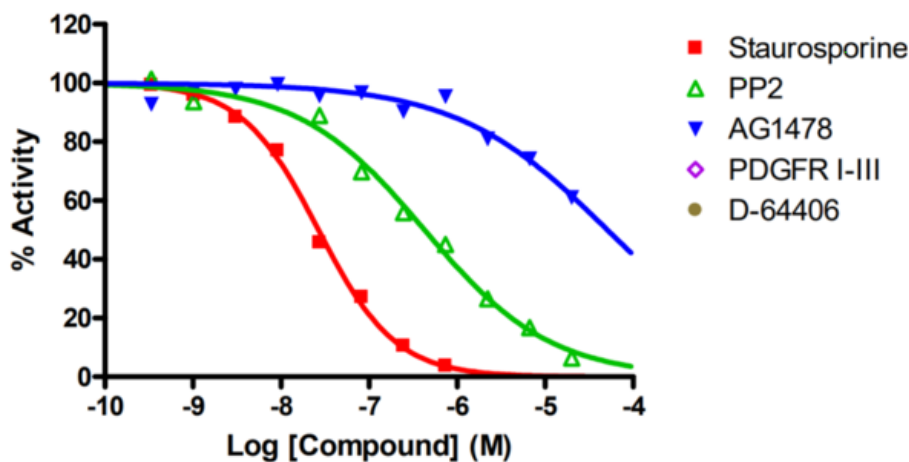
Target Overview

| | |
|------------|--|
| CAT# | BTK (E41K) |
| Enzyme | Human BTK (E41K) |
| HGNC | BTK (E41K) |
| Synonym | AGMX1, AT, ATK, BPK, PSCTK1 |
| 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 | <table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>26.7</td> </tr> <tr> <td>PP2</td> <td>428</td> </tr> <tr> <td>AG1478</td> <td>51,000</td> </tr> </tbody> </table> | Compounds | IC50 (nM) | Staurosporine | 26.7 | PP2 | 428 | AG1478 | 51,000 |
| Compounds | IC50 (nM) | | | | | | | | |
| Staurosporine | 26.7 | | | | | | | | |
| PP2 | 428 | | | | | | | | |
| AG1478 | 51,000 | | | | | | | | |
| Scientific Information | Q06187 (http://www.uniprot.org/uniprot/Q06187) | | | | | | | | |
| Screening Location | Malvern, USA | | | | | | | | |

Reference compound IC50 for BTK (E41K)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/btk_e41k_kin_malvern

