

TAK1 Kinase Assay Service

Activity-based TAK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

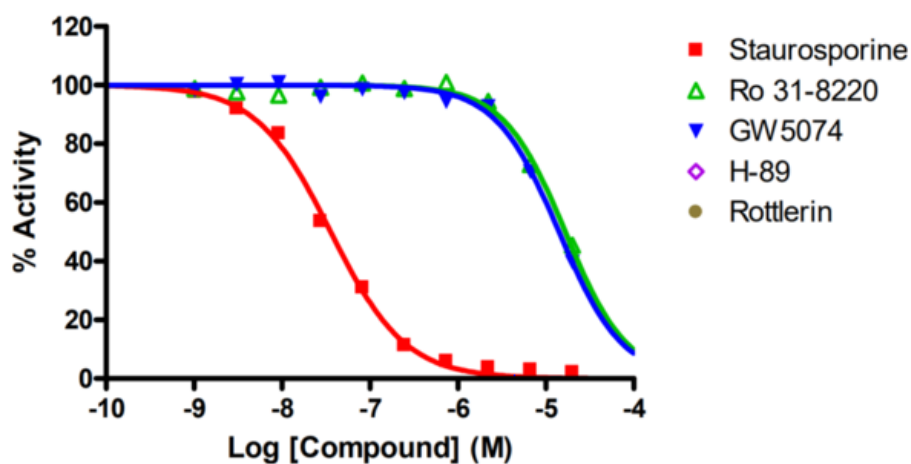
Target Overview

CAT#	TAK1
Enzyme	Human TAK1
HGNC	MAP3K7
Synonym	TGF1a, MEKK7
Substrates	Casein, 1 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>36</td> </tr> <tr> <td>Ro 31-8220</td> <td>16,700</td> </tr> <tr> <td>GW 5074</td> <td>14,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	36	Ro 31-8220	16,700	GW 5074	14,000
Compounds	IC50 (nM)								
Staurosporine	36								
Ro 31-8220	16,700								
GW 5074	14,000								
Scientific Information	O43318 (http://www.uniprot.org/uniprot/O43318)								
Screening Location	Malvern, USA								

Reference compound IC50 for TAK1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/tak1_kin_malvern

