

# ALK6/BMPR1B Kinase Assay Service

Activity-based ALK6/BMPR1B kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

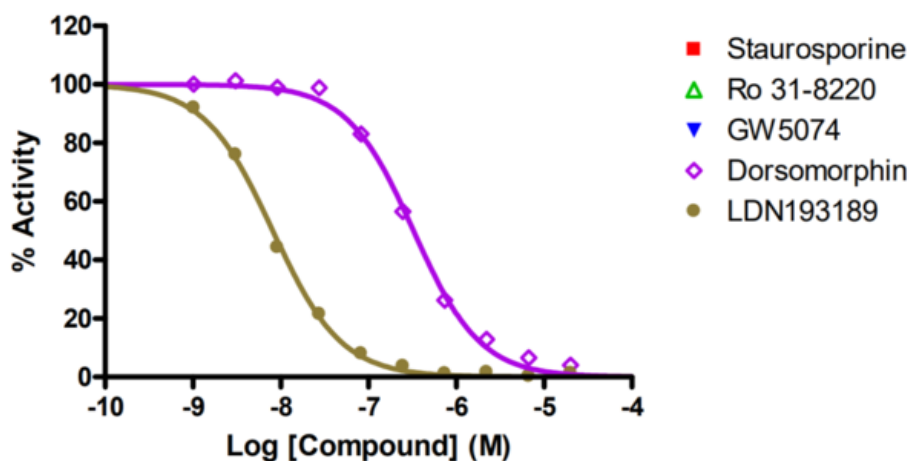
## Target Overview

CAT#	ALK6
Enzyme	Human ALK6
HGNC	BMPR1B
Synonym	ALK6, CDw293
Substrates	Casein, 20 μM ATP 10 μM

## Assay Properties

Reaction	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> 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>&gt;100,000</td> </tr> <tr> <td>Ro 31-8220</td> <td>&gt;100,000</td> </tr> <tr> <td>GW5074</td> <td>&gt;100,000</td> </tr> <tr> <td>Dorsomorphin</td> <td>340</td> </tr> <tr> <td>LDN193189</td> <td>9.4</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	>100,000	Ro 31-8220	>100,000	GW5074	>100,000	Dorsomorphin	340	LDN193189	9.4
Compounds	IC50 (nM)												
Staurosporine	>100,000												
Ro 31-8220	>100,000												
GW5074	>100,000												
Dorsomorphin	340												
LDN193189	9.4												
Scientific Information	<a href="http://www.uniprot.org/uniprot/O00238">O00238</a> ( <a href="http://www.uniprot.org/uniprot/O00238">http://www.uniprot.org/uniprot/O00238</a> )												
Screening Location	Malvern, USA												

## Reference compound IC50 for ALK6/BMPR1B



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/alk6\\_bmpr1b\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/alk6_bmpr1b_kin_malvern)

