

BPTF-[PHD-BRD] Bromodomain Assay Service (AlphaScreen)

BPTF-[PHD-BRD]-His [Reader Domain Assay](#) for compound screening and profiling using the AlphaScreen assay format to measure compound binding to the target.

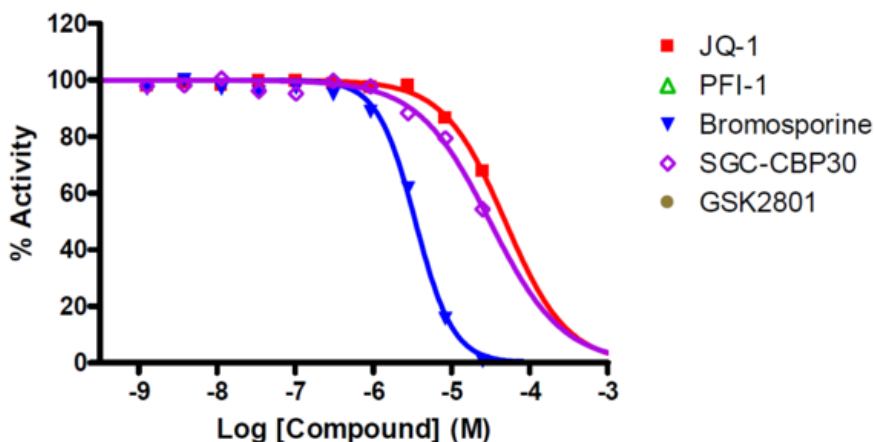
Target Overview

CAT#	BPTF-[PHD-BRD]-His
Enzyme	Human BPTF-[PHD-BRD]-His
Synonyms	Nucleosome-remodeling factor subunit BPTF, Bromodomain and PHD finger-containing transcription factor, Fetal Alz-50 clone 1 protein, Fetal Alzheimer antigen, FAC1, FALZ
Substrates	Histone H4 peptide (1-21) K5/8/12/16Ac

Assay Properties

Reaction	AlphaScreen Binding assay								
Readout	Fluorescence								
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>JQ-1</td> <td>50,000</td> </tr> <tr> <td>Bromosporine</td> <td>3,500</td> </tr> <tr> <td>SGC-CBP30</td> <td>31,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	JQ-1	50,000	Bromosporine	3,500	SGC-CBP30	31,000
Compounds	IC50 (nM)								
JQ-1	50,000								
Bromosporine	3,500								
SGC-CBP30	31,000								
Scientific Information	Q12830 (http://www.uniprot.org/uniprot/Q12830)								
Screening Location	Malvern, USA								
Further information	More information can be found on our website Reader Domain Assay Services .								

Reference Compound IC50 for BPTF-[PHD-BRD]



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/bptf-phd-brd-hisreader_alphascreen_malvern

