

# BRD2-2 Bromodomain Assay Service (AlphaScreen)

BRD2-2-His [Reader Domain Assay](#) for compound screening and profiling using the AlphaScreen assay format to measure compound binding to the target.

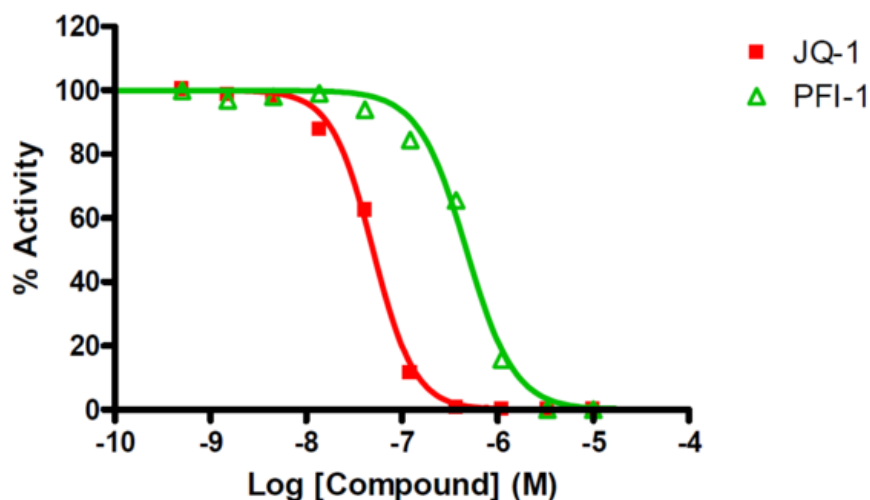
## Target Overview

CAT#	BRD2-2
Enzyme	Human BRD2-2
Synonyms	Bromodomain-containing protein 2 domain 2, O27.1.1, Really interesting new gene 3 protein, KIAA9001, RING3
Substrates	Histone H4 peptide (1-21) K5/8/12/16Ac

## Assay Properties

Reaction	AlphaScreen Binding assay
Readout	Fluorescence
Reference Compound IC50s	<b>Compounds IC50 (nM)</b> JQ-1      50 PFI-1     475
Scientific Information	<a href="http://www.uniprot.org/uniprot/P25440">P25440</a> ( <a href="http://www.uniprot.org/uniprot/P25440">http://www.uniprot.org/uniprot/P25440</a> )
Screening Location	Malvern, USA
Further information	More information can be found on our website <a href="#">Reader Domain Assay Services</a> .

## Reference Compound IC50 for BRD2-2



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/brd2-2reader\\_alphascreen\\_malvern](https://www.reactionbiology.com/datasheet/brd2-2reader_alphascreen_malvern)

