

BRD4 Assay Service (AlphaScreen)

BRD4 Full Length [Reader Domain Assay](#) for compound screening and profiling using the AlphaScreen assay format to measure compound binding to the target.

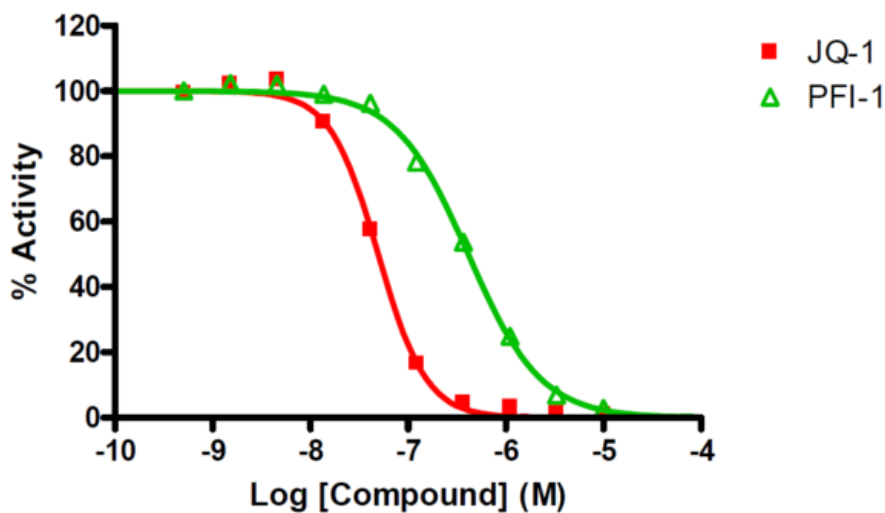
Target Overview

CAT#	BRD4 Full Length
Enzyme	Human BRD4 Full length
Synonyms	Bromodomain-containing protein 4 domain 1 and 2, Protein HUNK1
Substrates	Histone H4 peptide (1-21) K5/8/12/16Ac

Assay Properties

Reaction	AlphaScreen Binding assay
Readout	Fluorescence
Reference Compound IC50s	Compounds IC50 (nM) JQ-1 50 PFI-1 385
Scientific Information	O60885 (http://www.uniprot.org/uniprot/O60885)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

Reference Compound IC50 for BRD4 Full length



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brd4-full-lengthreader_alphascreen_malvern

