

BRD3 Bromodomain Assay Service (AlphaScreen)

BRD3-Tdndm-His [Reader Domain Assay](#) for compound screening and profiling using the AlphaScreen assay format to measure compound binding to the target.

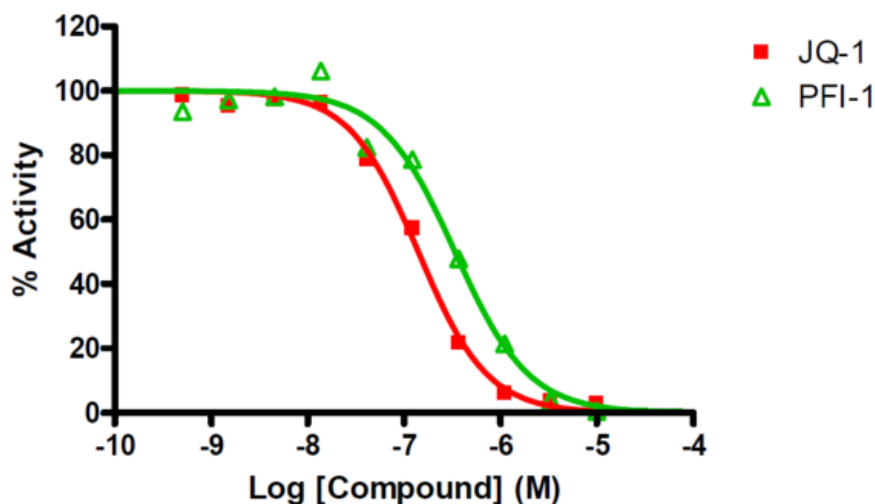
Target Overview

CAT#	BRD3-Tandem
Enzyme	Human BRD3-Tandem
Synonyms	Bromodomain-containing protein 3 domain 1 and domain 2, RING3-like protein, KIAA0043, RING3L
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-1 160 PFI-1 370
Scientific Information	Q15059 (http://www.uniprot.org/uniprot/Q15059)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

Reference Compound IC50 for BRD3-T



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brd3-tndmreader_alphascreen_malvern

