

SMARCA4 Bromodomain Assay Service (AlphaScreen)

SMARCA4-His [Reader Domain Assay](#) for compound screening and profiling using the AlphaScreen assay format to measure compound binding to the target.

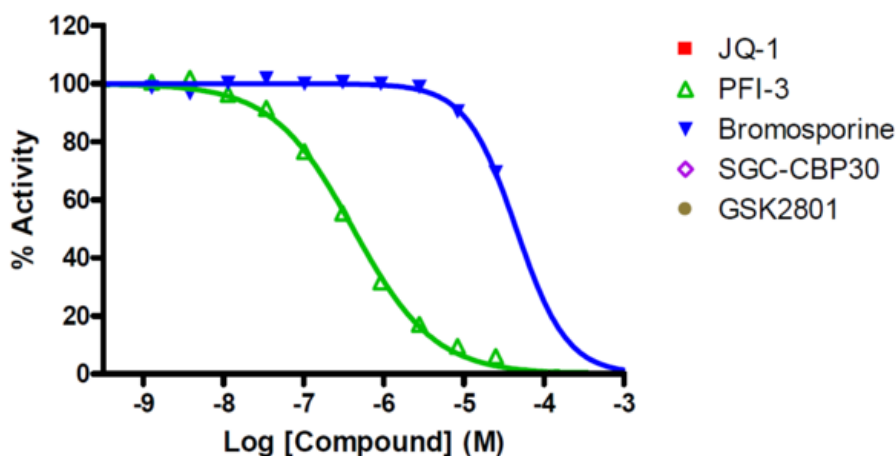
Target Overview

| | |
|------------|--|
| CAT# | SMARCA4-His |
| Enzyme | Human SMARCA4-His |
| Synonyms | Transcription activator BRG1, ATP-dependent helicase SMARCA4, BAF190A, BRG1, SNF2B, SNF2L4 |
| Substrates | Histone H3 peptide (1-30) K4/18/23/27Ac |

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>PFI-3</td> <td>420</td> </tr> <tr> <td>Bromosporine</td> <td>45,000</td> </tr> </tbody> </table> | Compounds | IC50 (nM) | PFI-3 | 420 | Bromosporine | 45,000 |
| Compounds | IC50 (nM) | | | | | | |
| PFI-3 | 420 | | | | | | |
| Bromosporine | 45,000 | | | | | | |
| Scientific Information | P51532 (http://www.uniprot.org/uniprot/P51532) | | | | | | |
| Screening Location | Malvern, USA | | | | | | |
| Further information | More information can be found on our website Reader Domain Assay Services . | | | | | | |

Reference Compound IC50 for SMARCA4



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/smarca4-hisreader_alphascreen_malvern

