

Bromosporine

Bromosporine is a synthetic small molecule pan-bromodomain inhibitor developed by the Structural Genomics Consortium for research. Used in research labs, and is shown to accelerate FRAP recovery of BRD4 and CREBBP in cells at 1uM concentration. Bromodomains are implicated in multiple cellular processes, including inflammatory gene expression, mitosis, and viral/host interaction.

Materials Provided

| | |
|---------------------------|---------------------------|
| Catalog Number: | 92-1239 |
| Quantity Supplied: | 1 vial(s), 10 mg per vial |

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

| | |
|--------------------------|---|
| Molecular Weight: | 404.44 g/mol |
| Source: | Synthetic |
| Purity: | ≥ 98% |
| Endotoxin Level: | N/A |
| Formulation: | N/A |
| Storage: | 2-8°C Please avoid multiple freeze/thaw cycles. |

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

| | |
|--------------------------------|----------|
| Stock Concentration: | 10 mM |
| Reconstitution Volume: | 2.473 mL |
| Reconstitution Solvent: | DMSO |

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com