

PFI-3

PFI-3 is a synthetic inhibitor of the SMARCA4 and PB1(5) bromodomains (KD=89 and 48nM, respectively). Also inhibits SMARCA2, and is otherwise selective over other bromodomain families. May play a role in SMARCA4 related cancers.

Materials Provided

| | |
|---------------------------|--------------------------|
| Catalog Number: | 92-1260 |
| Quantity Supplied: | 1 vial(s), 1 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: | 321.4 g/mol |
| Source: | Synthetic |
| Purity: | ≥ 98% |
| Endotoxin Level: | N/A |
| Formulation: | N/A |
| Storage: | -20°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: | 0.311 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