

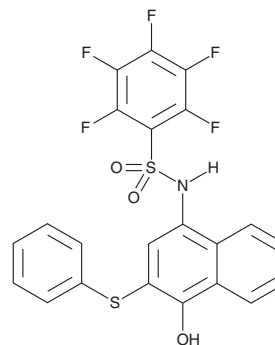
# PRODUCT INFORMATION



## MM-206

Item No. 22032

**CAS Registry No.:** 1809581-87-0  
**Formal Name:** 2,3,4,5,6-pentafluoro-N-[4-hydroxy-3-(phenylthio)-1-naphthalenyl]-benzenesulfonamide  
**MF:** C<sub>22</sub>H<sub>12</sub>F<sub>5</sub>NO<sub>3</sub>S<sub>2</sub>  
**FW:** 497.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 247 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

MM-206 is supplied as a crystalline solid. A stock solution may be made by dissolving the MM-206 in the solvent of choice, which should be purged with an inert gas. MM-206 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of MM-206 in these solvents is approximately 16 and 25 mg/ml, respectively.

MM-206 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MM-206 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MM-206 has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

MM-206 is an inhibitor of STAT3 (IC<sub>50</sub> = 1.16 μM in an SPR-based competitive assay).<sup>1</sup> It inhibits STAT3 phosphorylation induced by granulocyte colony-stimulating factor (G-CSF) in HL-60, Kasumi-1, and MOLM-13 acute myeloid leukemia (AML) cells (IC<sub>50</sub>s = 0.8, 1.9, and 1.1 μM, respectively). MM-206 induces apoptosis more potently in AML cell lines and patient-derived primary AML tumor cells than in acute lymphoblastic leukemia cells, which do not have upregulated STAT3 activity. It also slows disease progression and improves survival in an AML mouse xenograft model when administered at a dose of 30 mg/kg per day.

### Reference

1. Minus, M.B., Liu, W., Vohidov, F., *et al.* Rhodium(II) proximity-labeling identifies a novel target site on STAT3 for inhibitors with potent anti-leukemia activity. *Angew Chem. Int. Ed. Engl.* **54(44)**, 13085-13089 (2015).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/03/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM