

# PRODUCT INFORMATION



## STF-31

Item No. 11173

**CAS Registry No.:** 724741-75-7  
**Formal Name:** 4-[[[4-(1,1-dimethylethyl)phenyl]sulfonyl]amino]methyl]-N-3-pyridinyl-benzamide

**MF:** C<sub>23</sub>H<sub>25</sub>N<sub>3</sub>O<sub>3</sub>S

**FW:** 423.5

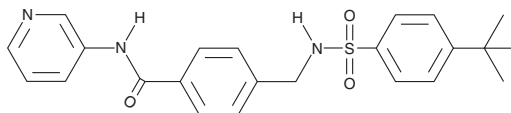
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 231, 264 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



### Laboratory Procedures

STF-31 is supplied as a crystalline solid. A stock solution may be made by dissolving the STF-31 in the solvent of choice. STF-31 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of STF-31 in these solvents is approximately 1, 20, and 25 mg/ml, respectively.

STF-31 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, STF-31 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. STF-31 has a solubility of approximately 0.2 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Glucose transporter 1 (Glut1) is an inducible carrier of pentoses and hexoses, including glucose. STF-31 is an inhibitor of Glut1 (IC<sub>50</sub> = ~1 μM) that blocks glucose uptake.<sup>1</sup> It induces necrosis in cancer cells that lack the von Hippel-Lindau tumor suppressor gene, which overexpress Glut1.<sup>1</sup> Although STF-31 binds Glut1, suggesting a direct effect, STF-31 also inhibits nicotinamide phosphoribosyltransferase, an enzyme that induces Glut1 expression.<sup>1-3</sup> STF-31 is also toxic to human pluripotent stem cells (hPSCs) and can be used to selectively eliminate hPSCs from mixed cultures.<sup>4</sup>

### References

1. Chan, D.A., Sutphin, P.D., Nguyen, P., *et al.* Targeting GLUT1 and the Warburg effect in renal cell carcinoma by chemical synthetic lethality. *Sci. Transl. Med.* **3(94)**, 1-9 (2011).
2. Adams, D.J., Ito, D., Rees, M.G., *et al.* NAMPT is the cellular target of STF-31-like small-molecule probes. *ACS Chem. Biol.* **9(10)**, 2247-2254 (2014).
3. Garten, A., Petzold, S., Körner, A., *et al.* Nampt: Linking NAD biology, metabolism, and cancer. *Trends Endocrinol. Metab.* **20(3)**, 130-138 (2009).
4. Boheler, K.R., Bhattacharya, S., Kropp, E.M., *et al.* A human pluripotent stem cell surface N-glycoproteome resource reveals markers, extracellular epitopes, and drug targets. *Stem Cell Reports* **3**, 185-203 (2014).

#### 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, 10/20/2016

#### 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