

# PRODUCT INFORMATION



## HA-130

Item No. 10498

**CAS Registry No.:** 1229652-21-4  
**Formal Name:** B-[3-[[4-[[3-[(4-fluorophenyl)methyl]-2,4-dioxo-5-thiazolidinylidene)methyl]phenoxy]methyl]phenyl]-boronic acid

**MF:** C<sub>24</sub>H<sub>19</sub>BFNO<sub>5</sub>S

**FW:** 463.3

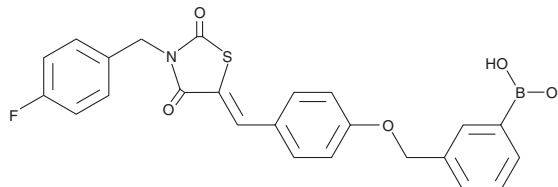
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 239, 348 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

HA-130 is supplied as a crystalline solid. A stock solution may be made by dissolving the HA-130 in the solvent of choice, which should be purged with an inert gas. HA-130 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of HA-130 in these solvents is approximately 0.5, 5, and 10 mg/ml, respectively.

### Description

HA-130 is a reversible inhibitor of autotaxin, completely blocking the hydrolysis of the substrate bis-pNPP with an IC<sub>50</sub> value of 28 nM.<sup>1</sup> It does not affect the activity of any proteasomal protease or related enzymes. HA-130 rapidly decreases plasma lysophosphatidic acid levels in mice when given intravenously (1 nM/g).<sup>1</sup> HA-130 has been used to investigate the role of autotaxin in cells and animals.<sup>2-4</sup>

### References

1. Albers, H. M. H. G., Dong, A., van Meeteren, L. A., *et al.* Boronic acid-based inhibitor of autotaxin reveals rapid turnover of LPA in the circulation. *Proc. Natl. Acad. Sci. USA* **107(16)**, 7257-7262 (2010).
2. Lai, S. L., Yao, W. L., Tsao, K. C., *et al.* Autotaxin/Lpar3 signaling regulates Kupffer's vesicle formation and left-right asymmetry in zebrafish. *Development* **139(23)**, 4439-4448 (2012).
3. Vázquez-Medina, J. P., Dodia, C., Weng, L., *et al.* The phospholipase A2 activity of peroxiredoxin 6 modulates NADPH oxidase 2 activation via lysophosphatidic acid receptor signaling in the pulmonary endothelium and alveolar macrophages. *FASEB J.* **30(8)**, 2885-2898 (2016).
4. Zhang, Y., Chen, Y. C. M., Krummel, M. F., *et al.* Autotaxin through lysophosphatidic acid stimulates polarization, motility, and transendothelial migration of naive T cells. *J. Immunol.* **189(8)**, 3914-3924 (2012).

#### 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, 10/04/2022

#### 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