

# PRODUCT INFORMATION



## HA-155

Item No. 11034

CAS Registry No.: 1229652-22-5

Formal Name: B-[4-[[4-[[3-[(4-fluorophenyl)methyl]-2,4-dioxo-5-thiazolidinylidene]methyl]phenoxy]methyl]phenyl]-boronic acid

Synonym: Autotaxin Inhibitor IV

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

FW: 463.3

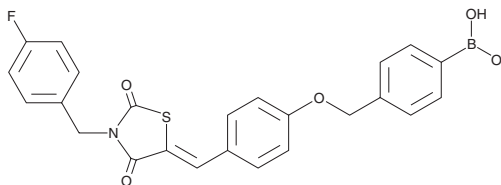
Purity: ≥95% (mixture of isomers)

UV/Vis.: λ<sub>max</sub>: 224, 347 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-155 is supplied as a crystalline solid. A stock solution may be made by dissolving the HA-155 in the solvent of choice, which should be purged with an inert gas. HA-155 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of HA-155 in these solvents is approximately 30 and 50 mg/ml, respectively.

HA-155 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Autotaxin converts lysophosphatidylcholine to lysophosphatidic acid (LPA), which can mediate changes in cell proliferation, angiogenesis, and cytokine secretion. HA-155 is a boronic acid-based compound that inhibits autotaxin (IC<sub>50</sub> = 5.7 nM) by selectively binding to its catalytic threonine.<sup>1,2</sup> It has been shown to dose-dependently block thrombin-induced LPA secretion in platelets.<sup>3</sup>

### References

1. Hausmann, J., Kamtekar, S., Christodoulou, E., *et al.* Structural basis of substrate discrimination and integrin binding by autotaxin. *Nat. Struct. Mol. Biol.* **18(2)**, 198-204 (2011).
2. Albers, H.M.H.G., Hendrickx, L.J.D., van Tol, R.J.P., *et al.* Structure-based design of novel boronic acid-based inhibitors of autotaxin. *J. Med. Chem.* **54(13)**, 4619-4626 (2011).
3. Fulkerson, Z., Wu, T., Sunkara, M., *et al.* Binding of autotaxin to integrins localizes lysophosphatidic acid production to platelets and mammalian cells. *J. Biol. Chem.* **286(40)**, 34654-34663 (2011).

#### 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, 11/14/2022

#### 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