

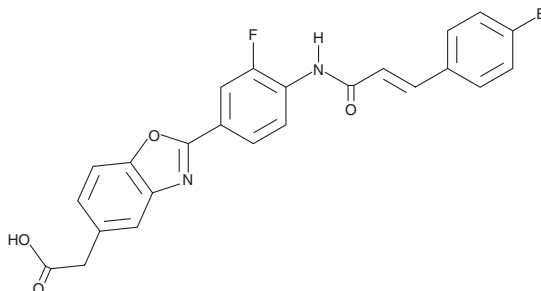
PRODUCT INFORMATION



OGT 2115

Item No. 17324

CAS Registry No.: 853929-59-6
Formal Name: 2-[4-[[3-(4-bromophenyl)-1-oxo-2-propen-1-yl]amino]-3-fluorophenyl]-5-benzoxazoleacetic acid
MF: C₂₄H₁₆BrFN₂O₄
FW: 495.3
Purity: ≥95%
Supplied as: A 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

OGT 2115 is supplied as a solid. A stock solution may be made by dissolving the OGT 2115 in the solvent of choice, which should be purged with an inert gas. OGT 2115 is soluble in DMSO.

Description

OGT 2115 is an inhibitor of heparanase (IC₅₀ = 0.4 μM), an enzyme that cleaves heparan sulfate into glucuronic acid (GlcUA) and N-acetylglucosamine (GlcNAc).^{1,2} OGT 2115 (10 μM) inhibits herpes simplex virus 2 (HSV-2) infection of VK2/E6E7 vaginal epithelial cells.³

References

1. Courtney, S.M., Hay, P.A., Buck, R.T., *et al.* Furanyl-1,3-thiazol-2-yl and benzoxazol-5-yl acetic acid derivatives: Novel classes of heparanase inhibitor. *Bioorg. Med. Chem. Lett.* **15(9)**, 2295-2299 (2005).
2. Nishimura, Y., Shitara, E., Adachi, H., *et al.* Flexible synthesis and biological activity of uronic acid-type gem-diamine 1-N-iminosugars: A new family of glycosidase inhibitors. *J. Org. Chem.* **65(1)**, 2-11 (2000).
3. Hopkins, J., Yadavalli, T., Agelidis, A.M., *et al.* Host enzymes heparanase and cathepsin L promote herpes simplex virus 2 release from cells. *J. Virol.* **92(23)**, e01179-e01218 (2018).

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, 04/06/2023

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