

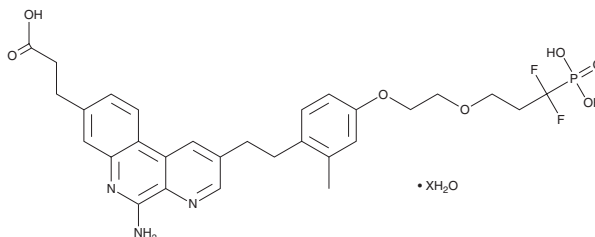
PRODUCT INFORMATION



LHC165 (hydrate)

Item No. 38776

Formal Name: 5-amino-2-[2-[4-[2-(3,3-difluoro-3-phosphonopropoxy)ethoxy]-2-methylphenyl]ethyl]-benzo[f][1,7]naphthyridine-8-propanoic acid, hydrate
MF: $C_{29}H_{32}F_2N_3O_7P \cdot XH_2O$
FW: 603.6
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 233 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

LHC165 (hydrate) is supplied as a solid. A stock solution may be made by dissolving the LHC165 (hydrate) in the solvent of choice, which should be purged with an inert gas. LHC165 (hydrate) is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

LHC165 is a toll-like receptor 7 (TLR7) agonist.¹

Reference

1. Zawit, M., Swami, U., Awada, H., *et al.* Current status of intralesional agents in treatment of malignant melanoma. *Ann. Transl. Med.* **9(12)**, 1038 (2021).

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, 07/24/2023

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