

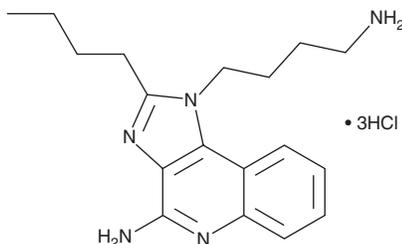
PRODUCT INFORMATION



AXC-715 (hydrochloride)

Item No. 44007

CAS Registry No.: 2479276-17-8
Formal Name: 4-amino-2-butyl-1H-imidazo[4,5-c]quinoline-1-butanamine, trihydrochloride
Synonym: T785
MF: C₁₈H₂₅N₅ • 3HCl
FW: 420.8
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

AXC-715 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the AXC-715 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AXC-715 (hydrochloride) is sparingly soluble (1-10 mg/ml) in the organic solvent DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AXC-715 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. AXC-715 (hydrochloride) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

AXC-715 is a dual agonist of toll-like receptor 7 (TLR7) and TLR8.¹ It activates NF-κB in a secreted alkaline phosphatase (SEAP) assay using TLR7- or TLR8-expressing HEK293 cells (EC₅₀s = 0.634 and 1.3 μM, respectively). AXC-715 has been linked to the anti-HER2 antibody trastuzumab as an antibody-drug conjugate (ADC). AXC-715 linked to trastuzumab reduces tumor volume in an HCC1954 breast cancer mouse xenograft model.

Reference

1. Ackerman, S.E., Pearson, C.I., Gregorio, J.D., *et al.* Immune-stimulating antibody conjugates elicit robust myeloid activation and durable antitumor immunity. *Nat. Cancer* **2**(1), 18-33 (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, 07/30/2025

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