

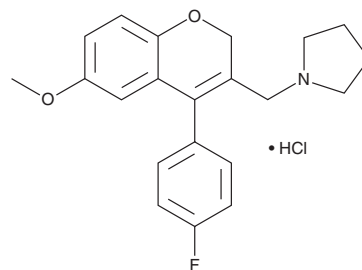
PRODUCT INFORMATION



AX-024 (hydrochloride)

Item No. 27333

CAS Registry No.: 1704801-24-0
Formal Name: 1-[[4-(4-fluorophenyl)-6-methoxy-2H-1-benzopyran-3-yl]methyl]-pyrrolidine, monohydrochloride
MF: C₂₁H₂₂FNO₂ • HCl
FW: 375.9
Purity: ≥98%
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

AX-024 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the AX-024 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AX-024 (hydrochloride) is soluble in the organic solvent DMSO.

Description

AX-024 is an inhibitor of the CD3ε-Nck interaction.¹ It binds to the SH3.1 N-terminal domain of Nck1, preventing its binding to the proline-rich sequence of the CD3ε subunit of T cells. AX-024 inhibits T cell receptor-triggered T cell activation (IC₅₀ = 1 nM for human T cells) and selectively decreases proliferation of T cells over B cells. It prevents CD4⁺ T cell and macrophage infiltration to the cerebellum and spinal cord and improves neurological impairments in a MOG₃₅₋₅₅ mouse model of experimental autoimmune encephalomyelitis (EAE) when administered at a dose of 10 mg/kg per day for ten days starting on the day of immunization. AX-024 (50 mg/kg) prevents the infiltration of eosinophils, macrophages, and neutrophils into the airway, as well as prevents an increase in IL-4 levels in bronchoalveolar lavage fluid (BALF), in an ovalbumin-sensitized mouse model of allergic airway disease. It does not prevent the memory T cell response to an immunodominant B8R poxvirus antigen peptide or to infection by ectromelia virus (mousepox).

Reference

1. Borroto, A., Reyes-Garau, D., Jiménez, M.A., *et al.* First-in-class inhibitor of the T cell receptor for the treatment of autoimmune diseases. *Sci. Transl. Med.* **8(370)**, 370ra184 (2016).

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, 12/08/2022

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