

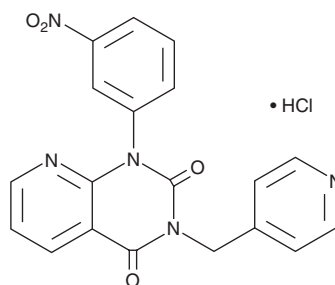
PRODUCT INFORMATION



RS 25344 (hydrochloride)

Item No. 37793

CAS Registry No.: 152815-28-6
Formal Name: 1-(3-nitrophenyl)-3-(4-pyridinylmethyl)-pyrido[2,3-d]pyrimidine-2,4(1H,3H)-dione
MF: C₁₉H₁₃N₅O₄ • HCl
FW: 411.8
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

RS 25344 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the RS 25344 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RS 25344 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of RS 25344 (hydrochloride) in these solvents is approximately 15 and 5 mg/ml, respectively.

Description

RS 25344 is an inhibitor of phosphodiesterase 4 (PDE4; IC₅₀ = 0.3 nM).¹ It is selective for PDE4 over PDE1, -2, and -3 (IC₅₀s = >100 μM for all). RS 25344 inhibits concanavalin A-induced IL-5 release and LPS-induced TNF-α release in isolated human peripheral blood mononuclear cells (PBMCs; EC₅₀s = 0.3 and 5.4 nM, respectively).²

References

1. Alvarez, R., Sette, C., Yang, D., *et al.* Activation and selective inhibition of a cyclic AMP-specific phosphodiesterase, PDE-4D3. *Mol. Pharmacol.* **48**(4), 616-622 (1995).
2. Draheim, R., Egerland, U., and Rundfeldt, C. Anti-inflammatory potential of the selective phosphodiesterase 4 inhibitor N-(3,5-dichloro-pyrid-4-yl)-[1-(4-fluorobenzyl)-5-hydroxy-indole-3-yl]-glyoxylic acid amide (AWD 12-281), in human cell preparations. *J. Pharmacol. Exp. Ther.* **308**(2), 555-563 (2004).

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, 06/16/2023

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