

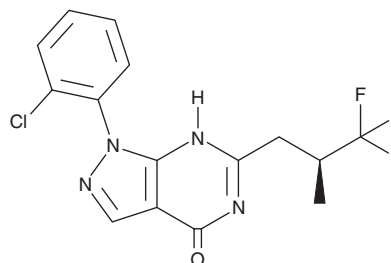
PRODUCT INFORMATION



(S)-BAY 73-6691

Item No. 44764

CAS Registry No.: 794568-91-5
Formal Name: 1-(2-chlorophenyl)-1,5-dihydro-6-
[(2S)-3,3,3-trifluoro-2-methylpropyl]-
4H-pyrazolo[3,4-d]pyrimidin-4-one
MF: C₁₅H₁₂ClF₃N₄O
FW: 356.7
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

(S)-BAY 73-6691 is supplied as a solid. A stock solution may be made by dissolving the (S)-BAY 73-6691 in the solvent of choice, which should be purged with an inert gas. (S)-BAY 73-6691 is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

(S)-BAY 73-6691 is an inhibitor of phosphodiesterase 9A (PDE9A; IC₅₀ = 88 nM) and the (S)-enantiomer of BAY 73-6691.¹

Reference

1. Wang, H., Luo, X., Ye, M., *et al.* Insight into binding of phosphodiesterase-9A selective inhibitors by crystal structures and mutagenesis. *J. Med. Chem.* **53**(4), 1726-1731 (2010).

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, 01/15/2026

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