

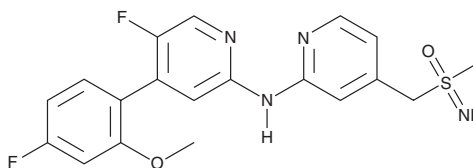
PRODUCT INFORMATION



(±)-BAY-1251152

Item No. 35094

CAS Registry No.: 1610358-53-6
Formal Name: 5-fluoro-4-(4-fluoro-2-methoxyphenyl)-N-[4-[(S-methylsulfonimidoyl)methyl]-2-pyridinyl]-2-pyridinamine
MF: C₁₉H₁₈F₂N₄O₂S
FW: 404.4
Purity: ≥98%
UV/Vis.: λ_{max}: 270, 329 nm
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

(±)-BAY-1251152 is supplied as a solid. A stock solution may be made by dissolving the (±)-BAY-1251152 in the solvent of choice, which should be purged with an inert gas. (±)-BAY-1251152 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (±)-BAY-1251152 in these solvents is approximately 12 mg/ml.

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 (±)-BAY-1251152 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of (±)-BAY-1251152 in PBS (pH 7.2) is approximately 0.16 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-BAY-1251152 is a racemic mixture of the cyclin-dependent kinase 9 (Cdk9) inhibitor (+)-BAY-1251152 and its isomer (-)-BAY-1251152.

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, 11/04/2022

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