

PRODUCT INFORMATION



BMS 345541 (trifluoroacetate salt)

Item No. 16667

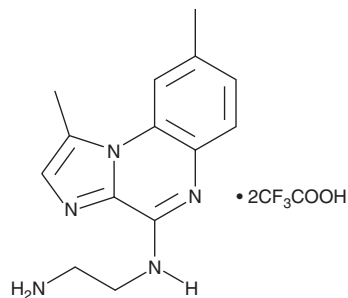
CAS Registry No.: 2320261-79-6
Formal Name: N¹-(1,8-dimethylimidazo[1,2-a]quinoxalin-4-yl)-1,2-ethanediamine, bis(2,2,2-trifluoroacetate)

Synonym: IKK Inhibitor III
MF: C₁₄H₁₇N₅ • 2CF₃COOH
FW: 483.4
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 226, 241, 261, 271, 291, 303, 316, 330 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BMS 345541 (trifluoroacetate salt) is supplied as a crystalline solid. Solutions of BMS 345541 (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of BMS 345541 (trifluoroacetate salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BMS 345541 is a cell permeable inhibitor of the IκB kinases IKKα and IKKβ (IC₅₀s = 4 and 0.3 μM).¹ It is without effect against a panel of other serine/threonine and tyrosine kinases at 100 μM.¹ BMS 345541 inhibits signaling through nuclear factor-κB (NF-κB) both in cells and *in vivo*, showing excellent pharmacokinetics in mice.¹ It blocks joint inflammation and damage in collagen-induced arthritis in mice and induces apoptosis in melanoma cells both *in vitro* and *in vivo*.^{2,3} BMS 345541 is used to explore novel roles for IKK phosphorylation and NF-κB signaling.^{4,5}

References

1. Burke, J.R., Pattoli, M.A., Gregor, K.R., *et al.* *J. Biol. Chem.* **278**(3), 1450-1456 (2003).
2. McIntyre, K.W., Shuster, D.J., Gillooly, K.M., *et al.* *Arthritis Rheum.* **48**(9), 2652-2659 (2003).
3. Yang, J., Amiri, K.I., Burke, J.R., *et al.* *Clin. Cancer Res.* **12**(3 Pt 1), 950-960 (2006).
4. Karim, Z.A., Zhang, J., Bangerjee, M., *et al.* *Blood* **121**(22), 4567-4574 (2013).
5. Wen, L., Zhu, M., Madigan, M.C., *et al.* *PLoS One* **9**(7), 1-12 (2014).

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, 02/12/2020

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