

PRODUCT INFORMATION



BTT3033

Item No. 42119

CAS Registry No.: 1259028-99-3
Formal Name: 1-(4-fluorophenyl)-N-methyl-N-[4-[[[(phenylamino)carbonyl]amino]phenyl]-1H-pyrazole-4-sulfonamide

MF: C₂₃H₂₀FN₅O₃S

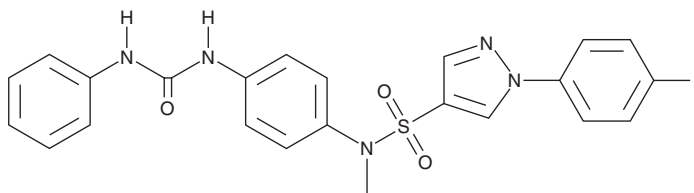
FW: 465.5

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

BTT3033 is supplied as a solid. A stock solution may be made by dissolving the BTT3033 in the solvent of choice, which should be purged with an inert gas. BTT3033 is soluble (≥10 mg/ml) in the organic solvent DMSO.

Description

BTT3033 is an inhibitor of integrin $\alpha 2\beta 1$ (EC_{50} = 130 nM in CHO cells expressing the human receptor).¹ It is 8-fold selective for integrin $\alpha 2\beta 1$ over integrin $\alpha 1\beta 1$, as well as laminin 332, 40 kDa fibronectin, 120 kDa fibronectin, and vitronectin at 130 nM. BTT3033 inhibits the adhesion of platelets in isolated mouse whole blood to collagen I-coated capillaries at a constant shear rate (EC_{50} = 6 μ M). *In vivo*, BTT3033 (10 mg/kg, p.o.) decreases lavage fluid leukocyte infiltration in a mouse model of platelet-activating factor-induced air pouch inflammation.² It decreases ear thickness in a mouse model of ovalbumin- and alum-induced mouse model of skin hypersensitivity when administered at a dose of 10 mg/kg.

References

1. Nissinen, L., Koivunen, J., Käpylä, J., *et al.* Novel $\alpha 2\beta 1$ integrin inhibitors reveal that integrin binding to collagen under shear stress conditions does not require receptor preactivation. *J. Biol. Chem.* **287**(53), 44694-44702 (2012).
2. Nissinen, L., Ojala, M., Langen, B., *et al.* Sulfonamide inhibitors of $\alpha 2\beta 1$ integrin reveal the essential role of collagen receptors in *in vivo* models of inflammation. *Pharmacol. Res. Perspect.* **3**(3), e00146 (2015).

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, 10/16/2024

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