

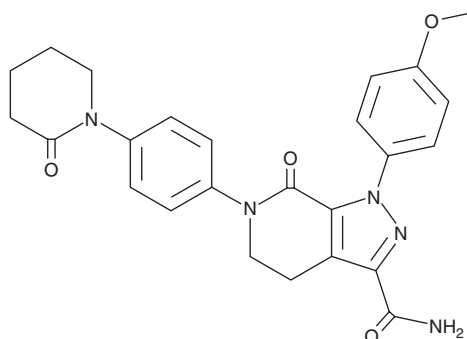
PRODUCT INFORMATION



Apixaban

Item No. 15427

CAS Registry No.: 503612-47-3
Formal Name: 4,5,6,7-tetrahydro-1-(4-methoxyphenyl)-7-oxo-6-[4-(2-oxo-1-piperidiny)phenyl]-1H-pyrazolo[3,4-c]pyridine-3-carboxamide
Synonym: BMS 562247-01
MF: C₂₅H₂₅N₅O₄
FW: 459.5
Purity: ≥98%
UV/Vis.: λ_{max}: 281 nm
Supplied as: A crystalline 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

Apixaban is supplied as a crystalline solid. A stock solution may be made by dissolving the apixaban in the solvent of choice, which should be purged with an inert gas. Apixaban is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of apixaban in these solvents is approximately 5 and 3 mg/ml, respectively.

Apixaban is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, apixaban should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Apixaban has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Apixaban is an orally bioavailable, selective inhibitor of both free and prothrombinase-bound factor Xa (K_s = 0.8 nM).¹ *In vivo*, apixaban has antithrombotic effects in a rabbit model of venous thrombosis. Apixaban (357.5 mg) also prevents thrombus formation without inducing adverse bleeding events in a porcine model of aortic heterotopic valve replacement.² Formulations containing apixaban have been used to prevent blood clot formation in patients with atrial fibrillation.

References

1. Turpie, A.G. Oral, direct factor Xa inhibitors in development for the prevention and treatment of thromboembolic diseases. *Arterioscler. Thromb. Vasc. Biol.* **27(6)**, 1238-1247 (2007).
2. Lester, P.A., Coleman, D.M., Diaz, J.A., et al. Apixaban versus warfarin for mechanical heart valve thromboprophylaxis in a swine aortic heterotopic valve model. *Arterioscler. Thromb. Vasc. Biol.* **37(5)**, 942-948 (2017).

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, 11/07/2022

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