

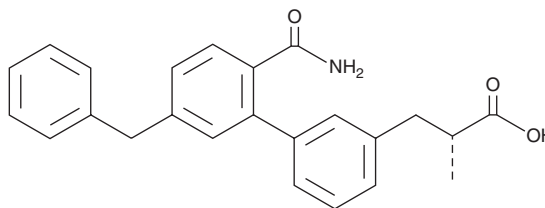
PRODUCT INFORMATION



AZD 2716

Item No. 35067

CAS Registry No.: 1845753-81-2
Formal Name: (αR)-2'-(aminocarbonyl)-α-methyl-5'-(phenylmethyl)-[1,1'-biphenyl]-3-propanoic acid
MF: C₂₄H₂₃NO₃
FW: 373.4
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

AZD 2716 is supplied as a solid. A stock solution may be made by dissolving the AZD 2716 in the solvent of choice, which should be purged with an inert gas. AZD 2716 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AZD 2716 in ethanol and DMF is approximately 16 mg/ml and approximately 10 mg/ml in DMSO.

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 AZD 2716 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of AZD 2716 in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 2716 is an inhibitor of secretory phospholipase A₂ (sPLA₂; IC₅₀s = 0.01, 0.04, and 0.4 μM for sPLA₂-IIa, -V, and -X, respectively).¹ It inhibits sPLA₂ activity in atherosclerotic plaque homogenates isolated from patients with coronary artery disease (IC₅₀ = 56 nM).

Reference

1. Giordanetto, F., Pettersen, D., Starke, I., *et al.* Discovery of AZD2716: A novel secreted phospholipase A₂ (sPLA₂) inhibitor for the treatment of coronary artery disease. *ACS Med. Chem. Lett.* **7**(10), 884-889 (2016).

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, 10/16/2025

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