

PRODUCT INFORMATION



AZD 7545

Item No. 19282

CAS Registry No.: 252017-04-2
Formal Name: 4-[[3-chloro-4-[[[(2R)-3,3,3-trifluoro-2-hydroxy-2-methyl-1-oxopropyl]amino]phenyl]sulfonyl]-N,N-dimethyl-benzamide

MF: C₁₉H₁₈ClF₃N₂O₅S
FW: 478.9

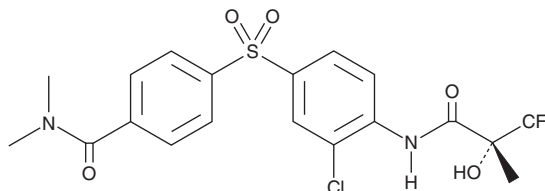
Purity: ≥98%

UV/Vis.: λ_{max}: 208, 273 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

AZD 7545 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 7545 in the solvent of choice, which should be purged with an inert gas. AZD 7545 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AZD 7545 in ethanol and DMF is approximately 20 mg/ml and approximately 50 mg/ml in DMSO.

AZD 7545 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 7545 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AZD 7545 has a solubility of approximately 0.02 mg/ml in a 1:40 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 7545 is an inhibitor of pyruvate dehydrogenase (PDH) kinases (PDKs), resulting in an increase in PDH activity with an EC₅₀ value of 5.2 nM for PDK2.^{1,2} It less potently inhibits PDK1 and PDK3 (IC₅₀s = 87 and 600 nM, respectively).³ AZD 7545 stimulates pyruvate oxidation in rat hepatocytes (EC₅₀ = 105 nM) and significantly elevates muscle PDH activity in obese Zucker rats.¹ It elevates postprandial glucose levels and improves the 24 hour glucose profile in obese, insulin-resistant rats, compared with lean counterparts.¹

References

1. Mayers, R.M., Butlin, R.J., Kilgour, E., *et al.* AZD7545, a novel inhibitor of pyruvate dehydrogenase kinase 2 (PDHK2), activates pyruvate dehydrogenase *in vivo* and improves blood glucose control in obese (*fa/fa*) Zucker rats. *Biochem. Soc. Trans.* **31(Pt 6)**, 1165-1167 (2003).
2. Morrell, J.A., Orme, J., Roche, T.E., *et al.* AZD7545 is a selective inhibitor of pyruvate dehydrogenase kinase 2. *Biochem. Soc. Trans.* **31(Pt 6)**, 1168-1170 (2016).
3. Kato, M., Li, J., Chuang, J.L., *et al.* Distinct structural mechanisms for inhibition of pyruvate dehydrogenase kinase isoforms by AZD7545, dichloroacetate, and radicicol. *Structure* **15(8)**, 992-1004 (2007).

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, 12/05/2022

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