

PRODUCT INFORMATION

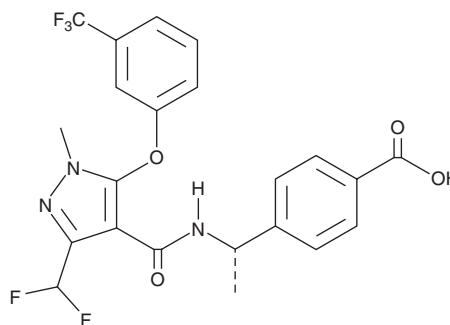


E7046

Item No. 19673

CAS Registry No.: 1369489-71-3
Formal Name: 4-[(1S)-1-[[[3-(difluoromethyl)-1-methyl-5-[3-(trifluoromethyl)phenoxy]-1H-pyrazol-4-yl]carbonyl]amino]ethyl]-benzoic acid

MF: C₂₂H₁₈F₅N₃O₄
FW: 483.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234 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

E7046 is supplied as a crystalline solid. A stock solution may be made by dissolving the E7046 in the solvent of choice, which should be purged with an inert gas. E7046 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of E7046 in these solvents is approximately 12, 14, and 2 mg/ml, respectively.

E7046 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, E7046 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. E7046 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

E7046 is an antagonist of the prostaglandin E₂ (PGE₂) receptor subtype EP₄ (IC₅₀ = 13.5 nM).¹ It inhibits PGE₂-induced activation of pro-tumor type 2 macrophages and prevents PGE₂-induced suppression of T cell proliferation in isolated human peripheral blood monocytes (PBMCs) when used at concentrations ranging from 0.1 to 3 μM. E7046 (5, 50, and 150 mg/kg) reduces tumor volume in Panc02 murine pancreatic and EMT6 murine breast cancer models.

Reference

1. Albu, D.I., Wang, Z., Huang, K.-C., *et al.* EP4 Antagonism by E7046 diminishes myeloid immunosuppression and synergizes with Treg-reducing IL-2-diphtheria toxin fusion protein in restoring anti-tumor immunity. *Oncoimmunology* **6**(8), e1338239 (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, 12/06/2022

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