

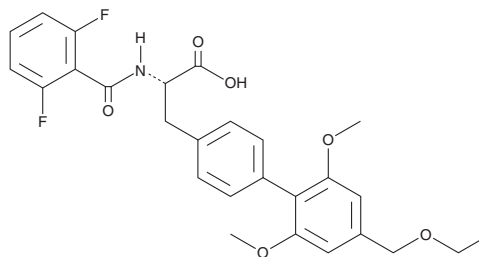
PRODUCT INFORMATION



Firategrast

Item No. 42893

CAS Registry No.: 402567-16-2
Formal Name: αS-[(2,6-difluorobenzoyl)amino]-4'-(ethoxymethyl)-2',6'-dimethoxy-[1,1'-biphenyl]-4-propanoic acid
Synonym: SB-683699
MF: C₂₇H₂₇F₂NO₆
FW: 499.5
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

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

Description

Firategrast is an antagonist of integrin α4β1, also known as very late antigen-4 (VLA-4), and integrin α4β7.¹ Maternal administration of firategrast (100 mg/kg) increases the number of donor hematopoietic stem cells (HSCs) in fetal liver in a mouse model of *in utero* allogeneic HSC transplant.

Reference

1. Kim, A.G., Vrecenak, J.D., Boelig, M.M., *et al.* Enhanced *in utero* allogeneic engraftment in mice after mobilizing fetal HSCs by α4β1/7 inhibition. *Blood* **128**(20), 2457-2461 (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, 02/12/2025

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