

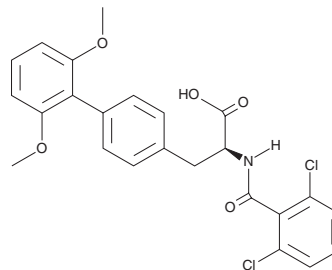
PRODUCT INFORMATION



TR-14035

Item No. 42726

CAS Registry No.: 232271-19-1
Formal Name: α S-[(2,6-dichlorobenzoyl)amino]-2',6'-dimethoxy-[1,1'-biphenyl]-4-propanoic acid
MF: C₂₄H₂₁Cl₂NO₅
FW: 474.3
Purity: $\geq 98\%$
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

TR-14035 is supplied as a solid. A stock solution may be made by dissolving the TR-14035 in the solvent of choice, which should be purged with an inert gas. TR-14035 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

TR-14035 is a dual $\alpha 4\beta 7$ and $\alpha 4\beta 1$ integrin antagonist.¹ It inhibits cell adhesion induced by the $\alpha 4\beta 7$ and $\alpha 4\beta 1$ integrin ligand fibronectin connecting segment 1 (FNCS1) in RPMI-8866 and Jurkat cells, which endogenously express $\alpha 4\beta 7$ and $\alpha 4\beta 1$, respectively (IC₅₀s = 7 and 87 nM, respectively).

Reference

1. Sircar, I., Gudmundsson, K.S., Martin, R., *et al.* Synthesis and SAR of N-benzoyl-L-biphenylalanine derivatives: Discovery of TR-14035, a dual $\alpha 4\beta 7/\alpha 4\beta 1$ integrin antagonist. *Bioorg. Med. Chem.* **10**(6), 2051-2066 (2002).

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, 02/18/2025

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