

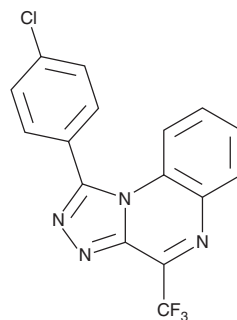
PRODUCT INFORMATION



R-8507

Item No. 22479

CAS Registry No.: 338773-13-0
Formal Name: 1-(4-chlorophenyl)-4-(trifluoromethyl)-[1,2,4]triazolo[4,3-a]quinoxaline
MF: C₁₆H₈ClF₃N₄
FW: 348.7
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 247 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

R-8507 is supplied as a crystalline solid. A stock solution may be made by dissolving the R-8507 in the solvent of choice, which should be purged with an inert gas. R-8507 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of R-8507 in these solvents is approximately 2 and 16 mg/ml, respectively.

R-8507 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, R-8507 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. R-8507 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

R-8507 is a small molecule antagonist of the TNF-α type 1 receptor (TNF-αRI).¹ It inhibits the expression of intercellular adhesion molecule-1 (ICAM-1) induced by TNF-α and IL-1β with EC₅₀ values of 2.45 and 3.79 μM, respectively, in an ELISA using A549 lung epithelial cells. It also disrupts the interaction of the TNF-αRI with receptor interacting protein 1 (RIP1) and TNF-αR-associated death domain protein (TRADD), preventing internalization of the receptor complex.

Reference

1. Gururaja, T.L., Yung, S., Ding, R., *et al.* A class of small molecules that inhibit TNFα-induced survival and death pathways via prevention of interactions between TNFαRI, TRADD, and RIP1. *Chemistry & Biology* **14**(10), 1105-1118 (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, 11/14/2022

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