

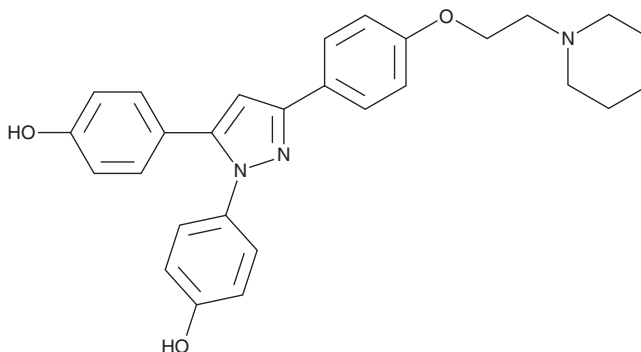
PRODUCT INFORMATION



TSI-13-57

Item No. 45498

CAS Registry No.: 2257427-45-3
Formal Name: 4,4'-[3-[4-[2-(1-piperidinyl)ethoxy]phenyl]-1H-pyrazole-1,5-diyl]bis-phenol
MF: C₂₈H₂₉N₃O₃
FW: 455.6
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

TSI-13-57 is supplied as a solid. A stock solution may be made by dissolving the TSI-13-57 in the solvent of choice, which should be purged with an inert gas. TSI-13-57 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

TSI-13-57 is an inhibitor of MyD88 adapter-like protein (Mal), also known as toll/IL-1 receptor domain-containing adaptor protein (TIRAP), homodimerization and of toll-like receptor (TLR) signaling.¹ It inhibits the secretion of TNF-α induced by agonists of TLR9, -7, -2, or -4 in RAW 264.7 macrophages (IC₅₀s = 2.7, 6.03, 7.37, and 6.38 μM, respectively). TSI-13-57 (100 and 200 mg/kg) reduces LPS-induced increases in TNF-α levels in mouse plasma.

Reference

1. Ippagunta, S. K., Pollock, J. A., Sharma, N., *et al.* Identification of toll-like receptor signaling inhibitors based on selective activation of hierarchically acting signaling proteins. *Sci. Signal* **11**(543), eaaq1077 (2018).

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, 04/03/2026

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