

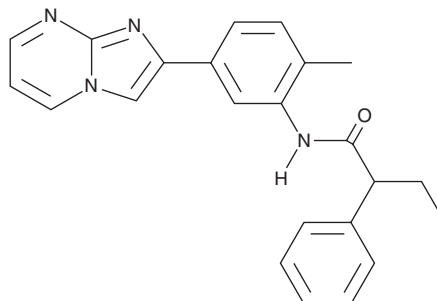
PRODUCT INFORMATION



PVP-037

Item No. 42073

CAS Registry No.: 862810-50-2
Formal Name: α -ethyl-N-(5-imidazo[1,2-a]pyrimidin-2-yl-2-methylphenyl)-benzeneacetamide
MF: $C_{23}H_{22}N_4O$
FW: 370.5
Purity: $\geq 95\%$
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

PVP-037 is supplied as a solid. A stock solution may be made by dissolving the PVP-037 in the solvent of choice, which should be purged with an inert gas. PVP-037 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PVP-037 is a toll-like receptor 7/8 (TLR7/8) agonist adjuvant.¹ It induces the production of cytokines and chemokines, including TNF, IFN- γ , IL-1 β , and IL-6, in primary human peripheral blood mononuclear cells (PBMCs) when used at concentrations greater than or equal to 1 μM . PVP-037, as an adjuvant to recombinant hemagglutinin protein (rHA), increases anti-rHA IgG titers in mice.

Reference

1. Soni, D., Borriello, F., Scott, D.A., et al. From hit to vial: Precision discovery and development of an imidazopyrimidine TLR7/8 agonist adjuvant formulation. Sci. Adv. 10(27), eadg3747 (2024).

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, 12/06/2024

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