

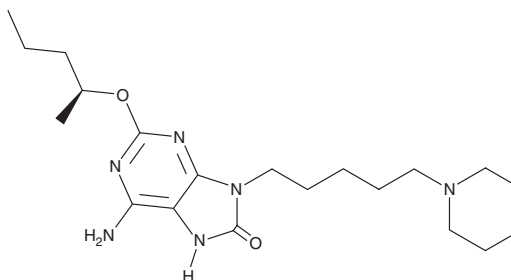
# PRODUCT INFORMATION



**GSK2245035**

Item No. 33171

**CAS Registry No.:** 1207629-49-9  
**Formal Name:** 6-amino-7,9-dihydro-2-[(1S)-1-methylbutoxy]-9-[5-(1-piperidinyl)pentyl]-8H-purin-8-one  
**MF:** C<sub>20</sub>H<sub>34</sub>N<sub>6</sub>O<sub>2</sub>  
**FW:** 390.5  
**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

GSK2245035 is supplied as a solid. A stock solution may be made by dissolving the GSK2245035 in the solvent of choice, which should be purged with an inert gas. GSK2245035 is soluble in the organic solvent DMSO. The solubility of GSK2245035 in DMSO is approximately 10 mM.

## Description

GSK2245035 is a toll-like receptor 7 (TLR7) agonist.<sup>1</sup> It selectively induces the production of IFN-α over TNF-α in isolated human peripheral blood mononuclear cells (PBMCs; EC<sub>50</sub>s = 0.63 and 316 nM, respectively), as well as over the production of TNF-α, IL-1β, IFN-γ, IL-10, and IL-12p70 in isolated human whole blood (EC<sub>50</sub>s = 63-10,000 nM). GSK2245035 (0.1-1,000 nM) reduces timothy grass- or house dust mite-induced IL-10 and IL-5 release in isolated human PBMCs. Intranasal administration of GSK2245035 (0.03-1 mg/kg) increases serum IP-10 levels, a marker of IFN-α induction, in mice. It also increases serum levels of IP-10 in cynomolgus monkeys.

## Reference

1. Biggadike, K., Ahmed, M., Ball, D.I., *et al.* Discovery of 6-amino-2-[[[(1S)-1-methylbutyl]oxy]-9-[5-(1-piperidinyl)pentyl]-7,9-dihydro-8H-purin-8-one (GSK2245035), a highly potent and selective intranasal toll-like receptor 7 agonist for the treatment of asthma. *J. Med. Chem.* **59**(5), 1711-1726 (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, 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](http://WWW.CAYMANCHEM.COM)