

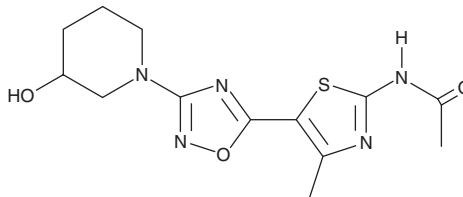
# PRODUCT INFORMATION



**TASP0415914**

Item No. 41145

**CAS Registry No.:** 1292300-75-4  
**Formal Name:** N-[5-[3-(3-hydroxy-1-piperidinyl)-1,2,4-oxadiazol-5-yl]-4-methyl-2-thiazolyl]-acetamide  
**MF:** C<sub>13</sub>H<sub>17</sub>N<sub>5</sub>O<sub>3</sub>S  
**FW:** 323.4  
**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

TASP0415914 is supplied as a solid. A stock solution may be made by dissolving the TASP0415914 in the solvent of choice, which should be purged with an inert gas. TASP0415914 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

TASP0415914 is a PI3K $\gamma$  inhibitor (IC<sub>50</sub> = 29 nM).<sup>1</sup> It is selective for PI3K $\gamma$  over Akt (IC<sub>50</sub> = 294 nM). *In vivo*, TASP0415914 (30 and 100 mg/kg) reduces ankle redness and swelling in a mouse model of collagen-induced arthritis.

## Reference

1. Oka, Y., Yabuuchi, T., Oi, T., *et al.* Discovery of N-{5-[3-(3-hydroxypiperidin-1-yl)-1,2,4-oxadiazol-5-yl]-4-methyl-1,3-thiazol-2-yl}acetamide (TASP0415914) as an orally potent phosphoinositide 3-kinase  $\gamma$  inhibitor for the treatment of inflammatory diseases. *Bioorg. Med. Chem.* **21**(24), 7578-7583 (2013).

### 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, 06/14/2024

## 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](http://WWW.CAYMANCHEM.COM)