

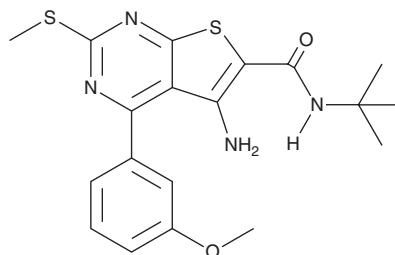
# PRODUCT INFORMATION



**ORG 41841**

Item No. 37724

**CAS Registry No.:** 301847-37-0  
**Formal Name:** 5-amino-N-(1,1-dimethylethyl)-4-(3-methoxyphenyl)-2-(methylthio)-thieno[2,3-d]pyrimidine-6-carboxamide  
**MF:** C<sub>19</sub>H<sub>22</sub>N<sub>4</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 402.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 242, 303 nm  
**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

ORG 41841 is supplied as a solid. A stock solution may be made by dissolving the ORG 41841 in the solvent of choice, which should be purged with an inert gas. ORG 41841 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 5 mg/ml. ORG 41841 is slightly soluble in ethanol and DMSO.

## Description

ORG 41841 is a dual luteinizing hormone (LH) and thyroid-stimulating hormone (TSH) receptor agonist.<sup>1</sup> It induces cAMP production in HEK293 cells expressing human LH or TSH receptors (EC<sub>50</sub>s = 0.3 and 6.5 μM, respectively).

## Reference

1. Moore, A., Jaeschke, H., Kleinau, G., *et al.* Evaluation of small-molecule modulators of the luteinizing hormone/choriogonadotropin and thyroid stimulating hormone receptors: Structure-activity relationships and selective binding patterns. *J. Med. Chem.* **49(13)**, 3888-3896 (2006).

### 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, 10/25/2022

## 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