

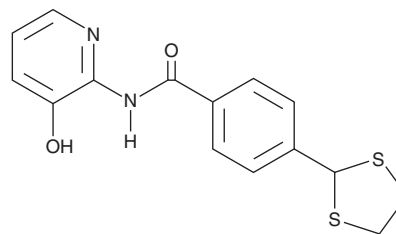
# PRODUCT INFORMATION



**E17241**

Item No. 41675

**CAS Registry No.:** 1060968-92-4  
**Formal Name:** 4-(1,3-dithiolan-2-yl)-N-(3-hydroxypyridin-2-yl)benzamide  
**Synonym:** 4-(1,3-Dithiolan-2-yl)-N-(3-hydroxypyridin-2-yl)benzamide  
**MF:** C<sub>15</sub>H<sub>14</sub>N<sub>2</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 318.4  
**Purity:** ≥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

E17241 is supplied as a solid. A stock solution may be made by dissolving the E17241 in the solvent of choice, which should be purged with an inert gas. E17241 is soluble (≥10 mg/ml) in DMSO.

## Description

E17241 is an inducer of the expression of ABCA1, the gene encoding ATP-binding cassette transporter 1; subfamily A (ABCA1).<sup>1</sup> It induces the expression of ABCA1 in a reporter assay using HepG2 cells (EC<sub>50</sub> = 280 nM). E17241 is also an agonist of peroxisome proliferator-activated receptors (PPARs).<sup>2</sup> It induces PPAR-mediated gene expression in reporter assays using HepG2 cells expressing PPAR $\gamma$ , PPAR $\alpha$ , or PPAR $\delta$  (EC<sub>50</sub>s = 290, 3,900, and 879 nM, respectively). It increases the protein levels of ACBA1 in RAW 264.7 macrophages but not in the presence of siRNA targeting mRNA encoding PKC $\zeta$ .<sup>1</sup> E17241 (0.4, 2, or 10  $\mu$ M) increases cholesterol efflux in RAW 264.7 cells. It reduces plasma levels of cholesterol, alanine transaminase (ALT), and aspartate aminotransferase (AST) and hepatic levels of cholesterol and triglycerides, as well as decreases aortic lesion area, in an ApoE<sup>-/-</sup> mouse model of atherosclerosis when administered at a dose of 25 mg/kg. E17241 (50 mg/kg per day) decreases plasma glucose levels and body weight in KKAY diabetic mice fed a high-fat and high-glucose (HFHG) diet.<sup>2</sup>

## References

- Xu, Y., Liu, C., Hang, X., *et al.* E17241 as a novel ABCA1 (ATP-binding cassette transporter A1) upregulator ameliorates atherosclerosis in mice. *Arterioscler. Thromb. Vasc. Biol.* **41(6)**, e284-e298 (2021).
- Sheng, R., Li, Y., Wu, Y., *et al.* A pan-PPAR agonist E17241 ameliorates hyperglycemia and diabetic dyslipidemia in KKAY mice via up-regulating ABCA1 in islet, liver, and white adipose tissue. *Biomed. Pharmacother.* **172**, 116200 (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, 10/16/2024

## 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