

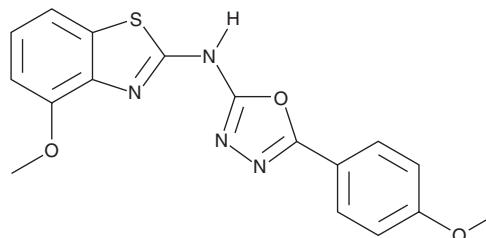
# PRODUCT INFORMATION



**N106**

Item No. 30952

**CAS Registry No.:** 862974-25-2  
**Formal Name:** 4-methoxy-N-[5-(4-methoxyphenyl)-1,3,4-oxadiazol-2-yl]-2-benzothiazolamine  
**MF:** C<sub>17</sub>H<sub>14</sub>N<sub>4</sub>O<sub>3</sub>S  
**FW:** 354.4  
**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

N106 is supplied as a solid. A stock solution may be made by dissolving the N106 in the solvent of choice, which should be purged with an inert gas. N106 is soluble in the organic solvent DMSO at a concentration of approximately 20 mg/ml.

## Description

N106 is an activator of sarcoplasmic/endoplasmic reticulum Ca<sup>2+</sup>-ATPase 2A (SERCA2A) SUMOylation.<sup>1</sup> It activates SUMO-activating enzyme E1, which increases SUMOylation and the activity of SERCA2A in isolated adult rat cardiomyocytes when used at concentrations ranging from 0.1 to 1 μM. N106 increases contractility in the same cells. *In vivo*, N106 (10 mg/kg) increases cardiac SERCA2A SUMOylation and contractility in a mouse model of pressure overload-induced heart failure.

## Reference

1. Kho, C., Lee, A., Jeong, D., *et al.* Small-molecule activation of SERCA2a SUMOylation for the treatment of heart failure. *Nat. Commun.* **6**, 7229 (2015).

### 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, 10/03/2022

## 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)