

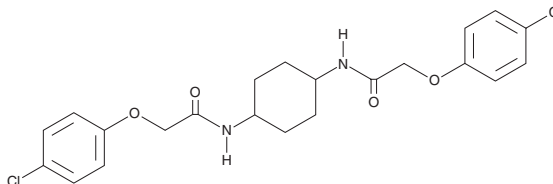
# PRODUCT INFORMATION



## ISRIB

Item No. 41879

CAS Registry No.: 548470-11-7  
Formal Name: N,N'-1,4-cyclohexanediylbis[2-(4-chlorophenoxy)-acetamide]  
MF: C<sub>22</sub>H<sub>24</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>4</sub>  
FW: 451.3  
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

ISRIB is supplied as a solid. A stock solution may be made by dissolving the ISRIB in the solvent of choice, which should be purged with an inert gas. ISRIB is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

### Description

ISRIB is a mixture of the diastereomers and integrated stress response inhibitors *cis*-ISRIB and *trans*-ISRIB (Item No. 16258).<sup>1</sup> *trans*-ISRIB inhibits gene transcription induced by activating transcription factor 4 (ATF4) more effectively than *cis*-ISRIB in a reporter assay using HEK293T cells (IC<sub>50</sub>s = 5 and 600 nM, respectively). *trans*-ISRIB decreases the latency to escape in the Morris water maze and increases the length of freezing time induced by contextual fear conditioning in mice.

### References

1. Sidrauski, C., Acosta-Alvear, D., Khoutorsky, A., *et al.* Pharmacological brake-release of mRNA translation enhances cognitive memory. *Elife* 2:e00498, (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, 11/05/2024

#### 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