

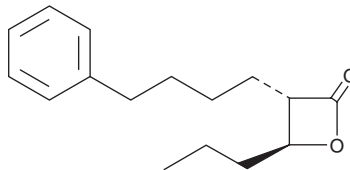
# PRODUCT INFORMATION



## (±)-*trans*-GK563

Item No. 25732

**CAS Registry No.:** 2351820-19-2  
**Formal Name:** (3R,4R)-*rel*-3-(4-phenylbutyl)-4-propyl-2-oxetanone  
**Synonym:** GK563  
**MF:** C<sub>16</sub>H<sub>22</sub>O<sub>2</sub>  
**FW:** 246.3  
**Purity:** ≥94%  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

(±)-*trans*-GK563 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. (±)-*trans*-GK563 is slightly soluble in chloroform, DMSO, and methanol.

### Description

*trans*-(±)-GK563 is a potent  $\beta$ -lactone inhibitor of group VIA (GVIA) calcium-independent phospholipase A<sub>2</sub> (GVIA iPLA<sub>2</sub>; IC<sub>50</sub> = 1 nM).<sup>1</sup> It is selective for GVIA iPLA<sub>2</sub> over GVIA cPLA<sub>2</sub> (IC<sub>50</sub> = 22  $\mu$ M). *trans*-(±)-GK563 (1 and 10  $\mu$ M) inhibits apoptosis induced by IL-1 $\beta$  and IFN- $\gamma$  in INS-1  $\beta$ -cells.

### Reference

1. Dedaki, C., Kokotou, M.G., Mouchlis, V.D., *et al.*  $\beta$ -Lactones: A novel class of Ca<sup>2+</sup>-independent phospholipase A<sub>2</sub> (group VIA iPLA<sub>2</sub>) inhibitors with the ability to inhibit  $\beta$ -cell apoptosis. *J. Med. Chem.* **62**(6), 2916-2927 (2019).

#### 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, 11/20/2019

#### 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