

# PRODUCT INFORMATION

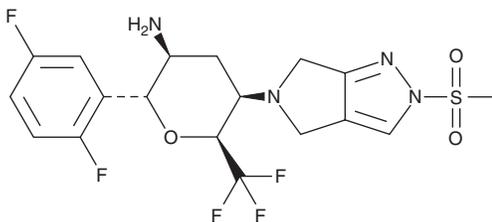


## Cofroglipitin

Item No. 45049

**CAS Registry No.:** 1844874-26-5  
**Formal Name:** 5-amino-2,6-anhydro-1,3,4,5-tetraoxo-6-C-(2,5-difluorophenyl)-3-[2,6-dihydro-2-(methylsulfonyl)pyrrolo[3,4-c]pyrazol-5(4H)-yl]-1,1,1-trifluoro-D-arabino-hexitol

**Synonyms:** HSK7653  
**MF:** C<sub>18</sub>H<sub>19</sub>F<sub>5</sub>N<sub>4</sub>O<sub>3</sub>S  
**FW:** 466.4  
**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

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

### Description

Cofroglipitin is a long-acting inhibitor of dipeptidyl peptidase 4 (DPP-4; IC<sub>50</sub> = 10.8 nM).<sup>1</sup> It is selective for DPP-4 over DPP-8 and DPP-9 (IC<sub>50</sub>s = >100 μM for both) and a panel of 127 receptors, ion channels, transporters, and enzymes at 10 μM. Cofroglipitin (10 mg/kg) decreases blood glucose levels, increases serum insulin levels, and inhibits serum DPP-4 activity in an oral glucose tolerance test using db/db mice.

### Reference

1. Gou, X., Dou, C., Tang, P., et al. Preclinical pharmacological profiles of cofroglipitin, a novel and bi-weekly DPP-4 inhibitor. *Front. Pharmacol.* **16**, 1702101 (2026).

**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, 02/06/2026

### 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