

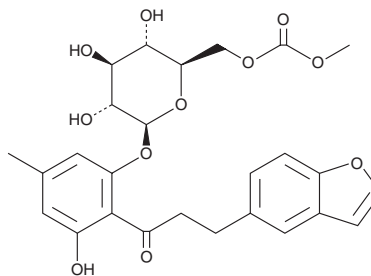
# PRODUCT INFORMATION



## T-1095

Item No. 37284

**CAS Registry No.:** 209746-59-8  
**Formal Name:** 3-(5-benzofuranyl)-1-[2-hydroxy-6-[[6-O-(methoxycarbonyl)-β-D-glucopyranosyl]oxy]-4-methylphenyl]-1-propanone  
**MF:** C<sub>26</sub>H<sub>28</sub>O<sub>11</sub>  
**FW:** 516.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

T-1095 is supplied as a crystalline solid. A stock solution may be made by dissolving the T-1095 in the solvent of choice, which should be purged with an inert gas. T-1095 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of T-1095 in these solvents is approximately 1, 2, and 5 mg/ml, respectively.

### Description

T-1095 is an inhibitor of sodium-glucose cotransporter 1 (SGLT1) and SGLT2 (IC<sub>50</sub>s = 22.8 and 2.3 μM, respectively).<sup>1</sup> Dietary administration of T-1095 (0.01% w/w) reduces blood glucose and HbA1c levels in a rat model of diabetes induced by streptozotocin (Item No. 13104). It suppresses the development of hyperglycemia and hypertriglyceridemia and reduces HbA1c levels in the KKAy mouse model of type 2 diabetes. T-1095 exacerbates post-infarction cardiac dysfunction in a rat model of left anterior descending coronary artery ligation-induced myocardial infarction.<sup>2</sup>

### References

1. Oku, A., Ueta, K., Ishihara, T., *et al.* T-1095, an inhibitor of renal Na<sup>+</sup>-glucose cotransporters, may provide a novel approach to treating diabetes. *Diabetes* **48(9)**, 1794-1800 (1999).
2. Bruno, J.G., Herman, T.S., Cano, V.L., *et al.* Selective cytotoxicity of 3-amino-L-tyrosine correlates with peroxidase activity. *In Vitro Cell. Dev. Biol. Anim.* **35(7)**, 376-382 (1999).

#### 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, 12/21/2022

#### 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