

PRODUCT INFORMATION



Telmisartan Acyl-β-D-Glucuronide (hydrate)

Item No. 28056

Formal Name: 1-[4'-[(1,4'-dimethyl-2'-propyl[2,6'-bi-1H-benzimidazol]-1'-yl)methyl][1,1'-biphenyl]-2-carboxylate] β-D-glucopyranuronic acid, hydrate

Synonym: Telmisartan Glucuronide

MF: C₃₉H₃₈N₄O₈ • XH₂O

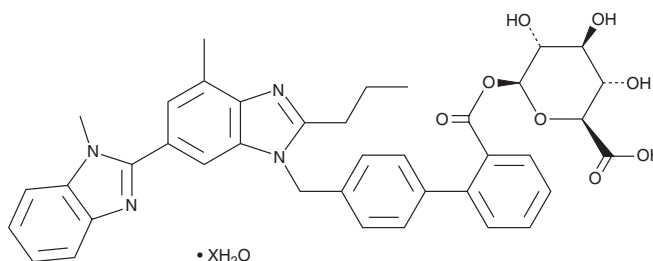
FW: 690.7

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

Telmisartan acyl-β-D-glucuronide (hydrate) is supplied as a solid. A stock solution may be made by dissolving the telmisartan acyl-β-D-glucuronide (hydrate) in the solvent of choice, which should be purged with an inert gas. Telmisartan acyl-β-D-glucuronide (hydrate) is soluble in methanol.

Description

Telmisartan acyl-β-D-glucuronide is a major metabolite of the angiotensin II receptor antagonist telmisartan (Item No. 11615).¹ It does not bind human serum albumin and is cleared from rat plasma with a clearance rate of 180 ml/min/kg following intravenous administration.

Reference

1. Ebner, T., Heinzl, G., Prox, A., *et al.* Disposition and chemical stability of telmisartan 1-O-acylglucuronide. *Drug Metab. Dispos.* **27**(10), 1143-1149 (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, 12/02/22

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