

# PRODUCT INFORMATION



## ZD 7155 (hydrochloride)

Item No. 36576

**CAS Registry No.:** 146709-78-6  
**Formal Name:** 5,7-diethyl-3,4-dihydro-1-[[2'-(2H-tetrazol-5-yl)[1,1'-biphenyl]-4-yl]methyl]-1,6-naphthyridin-2(1H)-one, monohydrochloride

**MF:** C<sub>26</sub>H<sub>26</sub>N<sub>6</sub>O • HCl

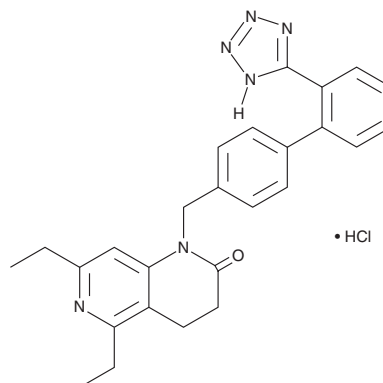
**FW:** 475.0

**Purity:** ≥98%

**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

ZD 7155 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ZD 7155 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ZD 7155 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

### Description

ZD 7155 is a nonpeptide antagonist of the angiotensin II type 1 (AT<sub>1</sub>) receptor (IC<sub>50</sub> = 3.3 nM).<sup>1</sup> It inhibits angiotensin II-induced pressor responses in conscious normotensive rats in a dose-dependent manner and decreases mean arterial pressure in conscious spontaneously hypertensive rats when administered at a dose of 1.082 μmol/kg.<sup>2</sup> ZD 7155 (10 mg/kg in the drinking water) improves motor nerve conduction velocity in the sciatic nerve in a rat model of streptozotocin-induced diabetic neuropathy.<sup>3</sup> It decreases disease severity and NF-κB and activator protein 1 (AP-1) activation in a mouse model of LPS-induced acute lung injury when administered at a dose of 10 mg/kg.<sup>4</sup>

### References

1. Zhang, H., Unal, H., Gati, C., *et al.* Structure of the Angiotensin receptor revealed by serial femtosecond crystallography. *Cell* **161**(4), 833-844 (2015).
2. Junggren, I.L., Zhao, X., Sun, X., *et al.* Comparative cardiovascular effects of the angiotensin II type 1 receptor antagonists ZD 7155 and losartan in the rat. *J. Pharm. Pharmacol.* **48**(8), 829-833 (1996).
3. Maxfield, E.K., Love, A., Cotter, M.A., *et al.* Nerve function and regeneration in diabetic rats: Effects of ZD-7155, an AT<sub>1</sub> receptor antagonist. *Am. J. Physiol.* **269**(3 Pt 1), E530-537 (1995).
4. Wang, F., Xia, Z.-F., Chen, X.-L., *et al.* Angiotensin II type-1 receptor antagonist attenuates LPS-induced acute lung injury. *Cytokine* **48**(3), 246-253 (2009).

#### 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, 12/20/2022

#### 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