

PRODUCT INFORMATION

DX600 (trifluoroacetate salt)

Item No. 22186

MF: $C_{141}H_{187}N_{35}O_{40}S_2 \cdot XCF_3COOH$

FW: 3,076.4

Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 221, 280 nm

Supplied as: A crystalline solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

Ac—Gly—Asp—Tyr—Ser—His—Cys—Ser—Pro—

Leu—Arg—Tyr—Tyr—Pro—Trp—Trp—Lys—Cys—

Thr—Tyr—Pro—Asp—Pro—Glu—Gly—Gly—Gly—NH₂

• XCF_3COOH

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

Laboratory Procedures

DX600 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the DX600 (trifluoroacetate salt) in the solvent of choice. DX600 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of DX600 (trifluoroacetate salt) is approximately 30 mg/ml in DMSO and DMF. DX600 is also slightly soluble in ethanol.

DX600 (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DX600 (trifluoroacetate salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DX600 (trifluoroacetate salt) has a solubility of approximately 0.13 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DX600 is a potent and selective peptide inhibitor of angiotensin-converting enzyme 2 (ACE2; K_i s = 2.8, 200, and 1,200 nM for recombinant, human, and murine ACE2, respectively).^{1,2} ACE2 is an enzyme that is cardioprotective and renoprotective.³ Inhibition of ACE2 with DX600 reduces *in vitro* vascular repair by CD34⁺ stem/progenitor cells and exacerbates myocardial injury and dysfunction *in vivo*.^{4,5}

References

- Huang, L., Sexton, D.J., Skogerson, K., *et al.* Novel peptide inhibitors of angiotensin-converting enzyme 2. *J. Biol. Chem.* **278**(18), 15532-15540 (2003).
- Joshi, S., Balasubramanian, N., Vasam, G., *et al.* Angiotensin converting enzyme versus angiotensin converting enzyme-2 selectivity of MLN-4760 and DX600 in human and murine bone marrow-derived cells. *Eur. J. Pharmacol.* **774**, 25-33 (2016).
- Prada, J. A. H., Ferreira, A. J., Katovich, M. J. *et al.* Structure-based identification of small-molecule angiotensin-converting enzyme 2 activators as novel antihypertensive agents. *Hypertension* **51**, 1312-1317 (2008).
- Singh, N., Joshi, S., Guo, L., *et al.* ACE2/Ang-(1-7)/Mas axis stimulates vascular repair-relevant functions of CD34⁺ cells. *Am. J. Physiol. Heart. Circ. Physiol.* **309**, H1697-H1707 (2015).
- Song, B., Zhang, Z.-Z., Zhong, J.-C., *et al.* Loss of angiotensin-converting enzyme 2 exacerbates myocardial injury via activation of the CTGF-fractalkine signaling pathway. *Circ. J.* **77**(12), 2997-3006 (2013).

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, 11/02/2022

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