

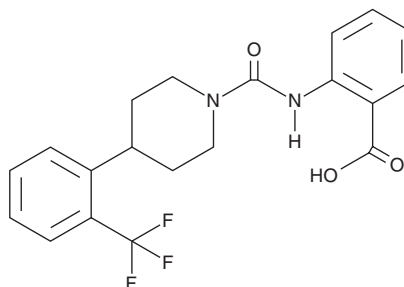
# PRODUCT INFORMATION



## A-1120

Item No. 30921

**CAS Registry No.:** 1152782-19-8  
**Formal Name:** 2-[[[4-[2-(trifluoromethyl)phenyl]-1-piperidiny]carbonyl]amino]-benzoic acid  
**MF:** C<sub>20</sub>H<sub>19</sub>F<sub>3</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 392.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 226, 252 nm  
**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

A-1120 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-1120 in the solvent of choice, which should be purged with an inert gas. A-1120 is soluble in ethanol and DMSO. The solubility of A-1120 in these solvents is approximately 25 and 100 mM, respectively.

### Description

A-1120 is a non-retinoid retinol binding protein 4 (RBP4) antagonist ( $K_i = 8.3$  nM for disruption of RBP4 binding to transthyretin).<sup>1</sup> It is selective for RBP4 over cellular retinol binding protein 1 (CRBP1;  $IC_{50}$ s = 0.09 and >30  $\mu$ M, respectively, in radioligand binding assays), as well as a panel of G protein-coupled receptors, kinases, and enzymes. A-1120 (30 mg/kg) reduces serum levels of RBP4 and retinol in mice. It also inhibits retinal bisretinoid accumulation in the *Abca4*<sup>-/-</sup> mouse model of enhanced retinal lipofuscinogenesis-induced macular degeneration.<sup>2</sup>

### References

1. Montani, A., Wang, Z., Conn, M., *et al.* Identification and characterization of a non-retinoid ligand for retinol-binding protein 4 which lowers serum retinol-binding protein 4 levels *in vivo*. *J. Biol. Chem.* **284**(12), 7673-7680 (2009).
2. Dobri, N., Qin, Q., Kong, J., *et al.* A1120, a nonretinoid RBP4 antagonist, inhibits formation of cytotoxic bisretinoids in the animal model of enhanced retinal lipofuscinogenesis. *Invest. Ophthalmol. Vis. Sci.* **54**(1), 85-95 (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, 11/01/2022

#### 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