

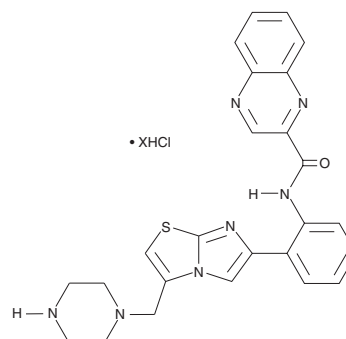
# PRODUCT INFORMATION



## SRT 1720 (hydrochloride)

Item No. 10011020

**CAS Registry No.:** 1001645-58-4  
**Formal Name:** N-[2-[3-(1-piperazinylmethyl)imidazo[2,1-b]thiazol-6-yl]phenyl]-2-quinolinecarboxamide, hydrochloride  
**Synonyms:** CAY10559, SRT 1933  
**MF:** C<sub>25</sub>H<sub>23</sub>N<sub>7</sub>OS • XHCl  
**FW:** 469.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 244 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

SRT 1720 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SRT 1720 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SRT 1720 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SRT 1720 (hydrochloride) in these solvents is approximately 10 and 1 mg/ml, respectively.

SRT 1720 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SRT 1720 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SRT 1720 (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Sirtuins (SIRT) represent a distinct class of trichostatin A-insensitive lysyl-deacetylases (class III HDACs). Human SIRT1 is the homolog of yeast silent information regulator 2 (Sir2) and has been shown to regulate the activity of the p53 tumor suppressor and inhibit apoptosis. Small molecule activators of SIRT1, such as resveratrol, extend lifespan in yeast and *C. elegans* in a manner that resembles caloric restriction. SRT 1720 is a selective small molecule activator of SIRT1 that is 1,000-fold more potent than resveratrol (EC<sub>1.5</sub> = 0.16 versus 46.2 μM, respectively).<sup>1</sup> In diet-induced obese and diabetic leptin-deficient *ob/ob* mice, oral administration of 100 mg/kg SRT 1720 once daily improves insulin sensitivity, lowers plasma glucose and increases mitochondrial capacity after one week of treatment.<sup>1</sup> In Zucker *fa/fa* rats, SRT 1720 improves whole-body glucose homeostasis and insulin sensitivity in adipose tissue, skeletal muscle, and liver.<sup>1</sup>

### Reference

1. Milne, J.C., Lambert, P.D., Schenk, S., *et al.* Small molecule activators of SIRT1 as therapeutics for the treatment of type 2 diabetes. *Nature* **450(7170)**, 712-716 (2007).

#### 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, 10/13/2022

#### 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