

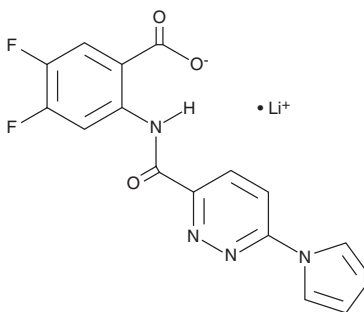
PRODUCT INFORMATION



SR 717 (lithium salt)

Item No. 31759

CAS Registry No.: 2375421-09-1
Formal Name: 4,5-difluoro-2-[[[6-(1H-imidazol-1-yl)-3-pyridazinyl]carbonyl]amino]-benzoic acid, monolithium salt
MF: C₁₅H₈F₂N₅O₃ • Li
FW: 351.2
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 290 nm
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

SR 717 (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the SR 717 (lithium salt) in the solvent of choice, which should be purged with an inert gas. SR 717 (lithium salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SR 717 (lithium salt) in these solvents is approximately 5 and 1 mg/ml, respectively.

SR 717 (lithium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SR 717 (lithium salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SR 717 (lithium salt) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SR 717 is an agonist of stimulator of interferon genes (STING; EC₅₀ = 2.1 μM in a reporter assay).¹ It is selective for STING over a panel of 468 kinases. SR 717 (30 mg/kg, i.p.) increases plasma IFN-β and IL-6 levels in mice. It also increases the intratumor frequency of CD69⁺CD8⁺ T cells, reduces tumor growth, and increases survival in a B16/F10 murine melanoma model.

Reference

1. Chin, E.N., Yu, C., Vartabedian, V.F., *et al.* Antitumor activity of a systemic STING-activating non-nucleotide cGAMP mimetic. *Science* **369**(6506), 993-999 (2020).

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, 10/19/2022

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