

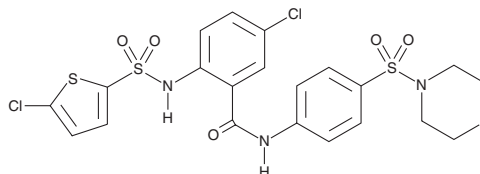
# PRODUCT INFORMATION



## Ataciguat

Item No. 16371

**CAS Registry No.:** 254877-67-3  
**Formal Name:** 5-chloro-2-[[[(5-chloro-2-thienyl)sulfonyl]amino]-N-[4-(4-morpholinylsulfonyl)phenyl]-benzamide  
**Synonym:** HMR 1766  
**MF:** C<sub>21</sub>H<sub>19</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>6</sub>S<sub>3</sub>  
**FW:** 576.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 241, 270 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

Ataciguat is supplied as a crystalline solid. A stock solution may be made by dissolving the ataciguat in the solvent of choice, which should be purged with an inert gas. Ataciguat is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 1 mg/ml.

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

### Description

Soluble guanylate cyclase (sGC) is the primary cellular receptor for NO. NO binds and activates a heme group in sGC, initiating the conversion of GTP to the second messenger cGMP. cGMP subsequently mediates a number of signaling cascades leading to vasorelaxation and inhibiting smooth muscle proliferation, leukocyte recruitment, and platelet aggregation. Oxidation of the heme results in its dissociation from sGC and an impairment of NO signaling, which has been linked to hypertension, hyperlipidemia, cardiovascular disease, and diabetes.<sup>1</sup> Ataciguat is an anthranilic acid derivative that activates the oxidized (heme-free) form of sGC (EC<sub>50</sub> = 0.5-10 μM in purified bovine lung or crude human corpus cavernosum isolates) by binding to the heme pocket and mimicking its function.<sup>1-3</sup> It has been shown to increase cGMP levels in cultured rat aorta smooth muscle cells and to induce vasorelaxation of isolated rat aorta, porcine coronary arteries, and human corpus cavernosum (EC<sub>50</sub> = 1-10 μM).<sup>2</sup>

### References

1. Stasch, J.P., Pacher, P., and Evgenov, O.V. Soluble guanylate cyclase as an emerging therapeutic target in cardiopulmonary disease. *Circulation* **123**(20), 2263-2273 (2011).
2. Schindler, U., Strobel, H., Schönafinger, K., et al. Biochemistry and pharmacology of novel anthranilic acid derivatives activating heme-oxidized soluble guanylyl cyclase. *Mol. Pharmacol.* **69**(4), 1260-1268 (2006).
3. Zhou, Z., Pyriochou, A., Kotanidou, A., et al. Soluble guanylyl cyclase activation by HMR-1766 (ataciguat) in cells exposed to oxidative stress. *Am. J. Physiol. Heart Circ. Physiol.* **295**(4), H1763-H1771 (2008).

#### 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, 02/04/2025

#### 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