

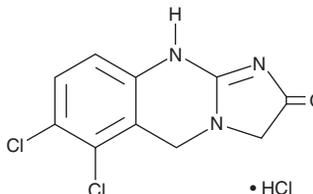
PRODUCT INFORMATION



Anagrelide (hydrochloride)

Item No. 21411

CAS Registry No.: 58579-51-4
Formal Name: 6,7-dichloro-1,5-dihydro-
imidazo[2,1-b]quinazolin-2(3H)-one,
monohydrochloride
Synonyms: BL 4162A, BMY 26538-01
MF: C₁₀H₇Cl₂N₃O • HCl
FW: 292.6
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 257 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

Anagrelide (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the anagrelide (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Anagrelide (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of anagrelide (hydrochloride) in these solvents is approximately 2 and 0.25 mg/ml, respectively.

Anagrelide (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, anagrelide (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Anagrelide (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

Anagrelide is an inhibitor of phosphodiesterase 3 (PDE3; IC₅₀ = 36 nM for the human platelet enzyme).¹ It inhibits thrombopoietin-induced megakaryocytopoiesis of isolated human CD34⁺ progenitor cells (IC₅₀ = 26 nM).² Anagrelide reduces platelet aggregation induced by ADP, collagen, thrombin, or arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) in isolated rabbit platelet rich-plasma (EC₅₀s = 0.31, 0.08, 0.18, and 0.1 μg/ml, respectively).³ It inhibits platelet thrombus formation in a dog model of electrically induced carotid artery thrombosis when administered at doses ranging from 0.5 to 5 mg/kg.⁴ Formulations containing anagrelide have been used in the treatment of thrombocytopenia.

References

1. Wang, G., Franklin, R., Hong, Y., *et al.* Comparison of the biological activities of anagrelide and its major metabolites in haematopoietic cell cultures. *Br. J. Pharmacol.* **146(3)**, 324-332 (2005).
2. Hong, Y., Wang, G., Gutierrez del Arroyo, A., *et al.* Comparison between anagrelide and hydroxycarbamide in their activities against haematopoietic progenitor cell growth and differentiation: Selectivity of anagrelide for the megakaryocytic lineage. *Leukemia* **20(6)**, 1117-1122 (2006).
3. Fleming, J.S. and Buyniski, J.P. A potent new inhibitor of platelet aggregation and experimental thrombosis, anagrelide (BL-4162A). *Thromb. Res.* **15(3-4)**, 373-388 (1979).
4. Fleming, J.S. and Buyniski, J.P. Anagrelide. *New drugs annual: Cardiovascular drugs*. Scriabine, A., editor, Raven Press (1983).

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, 01/04/2023

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