

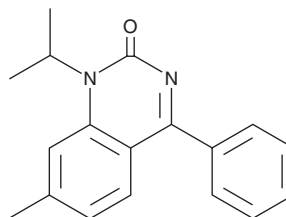
PRODUCT INFORMATION



Proquazone

Item No. 45454

CAS Registry No.: 22760-18-5
Formal Name: 7-methyl-1-(1-methylethyl)-4-phenyl-2(1H)-quinazolinone
Synonym: RU 43-715
MF: C₁₈H₁₈N₂O
FW: 278.4
Purity: ≥98%
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

Proquazone is supplied as a solid. A stock solution may be made by dissolving the proquazone in the solvent of choice, which should be purged with an inert gas. Proquazone is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Proquazone is a non-steroidal anti-inflammatory drug (NSAID).^{1,2} It reduces carrageenan-induced edema in rats (ED₅₀ = 5 mg/kg, p.o.) and bradykinin-induced bronchoconstriction in guinea pigs (ED₅₀ = 0.01 mg/kg, i.v.).¹ It also inhibits yeast-induced pyresis in rats.² Formulations containing proquazone have been used in the treatment of rheumatoid arthritis, osteoarthritis, and ankylosing spondylitis.

References

1. Coombs, R.V., Danna, R.P., Denzer, M., *et al.* Synthesis and antiinflammatory activity of 1-alkyl-4-aryl-2(1H)-quinazolinones and quinazolinethiones. *J. Med. Chem.* **16**(11), 1237-1245 (1973).
2. Ott, H. 2 (1H)-quinazolinones as novel non-acidic anti-inflammatory agents. *Scand. J. Rheumatol. Suppl.* **21**, 5-7 (1978).

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, 04/23/2026

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