

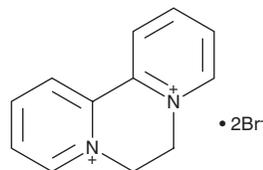
PRODUCT INFORMATION



Diquat (bromide)

Item No. 26251

CAS Registry No.: 85-00-7
Formal Name: 6,7-dihydro-dipyrido[1,2-a:2',1'-c]pyrazinediium, dibromide
MF: $C_{12}H_{12}N_2 \cdot 2Br$
FW: 344.0
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 312 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Diquat (bromide) is supplied as a crystalline solid. A stock solution may be made by dissolving the diquat (bromide) in the solvent of choice. Diquat (bromide) is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 1 mg/ml. Diquat (bromide) is also slightly soluble in ethanol.

Description

Diquat is a bipyridylum herbicide that, as a stable radical, acts as an electron acceptor in the electron transport chain and is reoxidized by molecular oxygen, producing reactive oxygen species (ROS).¹ Diquat administration has been used to induce oxidative stress *in vitro* and in animal models.²⁻⁴ It increases the production of ROS and decreases the mitochondrial membrane potential in mitochondria isolated from porcine intestine leading to mitophagy when administered at a dose of 100 mg/kg.² Formulations containing diquat have been used in agriculture for crop desiccation and defoliation.

References

1. Moreland, D.E. Mechanisms of action of herbicides. *Ann. Rev. Plant Physiol.* **31**, 597-638 (1980).
2. Cao, S., Wu, H., Wang, C., *et al.* Diquat-induced oxidative stress increases intestinal permeability, impairs mitochondrial function, and triggers mitophagy in piglets. *J. Anim. Sci.* **96(5)**, 1795-1805 (2018).
3. Singh, P., Hanson, P.S., and Morris, C.M. Sirtuin-2 protects neural cells from oxidative stress and is elevated in neurodegeneration. *Parkinsons Dis.* **2017:2643587**, (2017).
4. Tomášek, O., Gabrielová, B., Kačer, P., *et al.* Opposing effects of oxidative challenge and carotenoids on antioxidant status and condition-dependent sexual signalling. *Sci. Rep.* **6:23546**, (2016).

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, 10/27/2022

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