

# PRODUCT INFORMATION



## 360A (iodide)

Item No. 42501

CAS Registry No.: 737763-37-0

Formal Name: 3,3'-[2,6-pyridinediylbis(carbonylimino)]bis[1-methyl-quinolinium, diiodide

MF:  $C_{27}H_{23}N_5O_2 \cdot 2I$

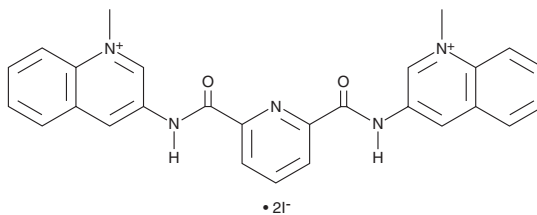
FW: 703.3

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^\circ\text{C}$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

360A (iodide) is supplied as a solid. A stock solution may be made by dissolving the 360A (iodide) in the solvent of choice, which should be purged with an inert gas. 360A (iodide) is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

360A is an inhibitor of telomerase ( $IC_{50} = 0.3 \mu\text{M}$ ).<sup>1</sup> It decreases the viability of in T98G glioma fibroblasts, CB193 astrocytoma and U-118 MG glioma cells, and primary human astrocytes ( $IC_{50}$ s = 4.8, 3.9, 8.4, and 17.4  $\mu\text{M}$ , respectively). 360A (5  $\mu\text{M}$ ) induces sister chromatid fusions and telomere loss in HeLa cervical cancer cells.<sup>2</sup> It induces cell cycle arrest at the S phase and increases  $\gamma$  histone H2AX ( $\gamma\text{H2AX}$ ) foci, a marker of DNA damage, in GM03657 B cell lymphocytes.

### References

1. Pennarun, G., Granotier, C., Gauthier, L.R., *et al.* Apoptosis related to telomere instability and cell cycle alterations in human glioma cells treated by new highly selective G-quadruplex ligands. *Oncogene* **24**(18), 2917-2928 (2005).
2. Pennarun, G., Granotier, C., Hoffschir, F., *et al.* Role of ATM in the telomere response to the G-quadruplex ligand 360A. *Nucleic Acids Res.* **36**(5), 1741-1754 (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, 02/12/2025

#### 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