# **PRODUCT** INFORMATION



GJ-103 (sodium salt)

Item No. 22482

CAS Registry No.:	1459687-96-7	
Formal Name:	2-[[4-(3-methoxyphenyl)-5-(2-	N
	pyridinyl)-4H-1,2,4-triazol-3-yl]	
	thio]-acetic acid, monosodium salt	
MF:	$C_{16}H_{13}N_4O_3S \bullet Na$	
FW:	364.4	N
Purity:	≥98%	
UV/Vis.:	λ <sub>max</sub> : 284 nm	·o
Supplied as:	A crystalline solid	S L
Storage:	-20°C	• Na+ 🛛 🗸
Stability:	≥4 years	

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

# Laboratory Procedures

GJ-103 (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the GJ-103 (sodium salt) in the solvent of choice, which should be purged with an inert gas. GJ-103 (sodium salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GJ-103 (sodium salt) in these solvents is approximately 30 mg/ml.

GJ-103 (sodium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GJ-103 (sodium salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GJ-103 (sodium salt) has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

# Description

GJ-103 is an inducer of stop codon read-through.<sup>1</sup> It restores the kinase activity of ataxia-telangiectasia mutated kinase (ATM) in patient-derived ataxia telangiectasia (A-T) lymphoblastoid cell lines containing homozygous TAG, TGA, or TAA nonsense mutations in ATM when used at concentrations of 10 and 30  $\mu$ M.

# Reference

1. Du, L., Jung, M.E., Damoiseaux, R., et al. A new series of small molecular weight compounds induce read through of all three types of nonsense mutations in the ATM gene. Mol. Ther. 21(9), 1653-1660 (2013).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

## SAFFTY 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

uyer 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, 11/22/2022

# 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