

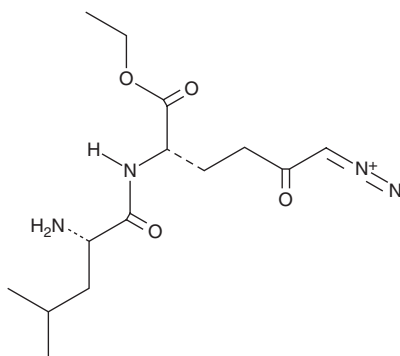
PRODUCT INFORMATION



JHU083

Item No. 40642

CAS Registry No.: 1998725-11-3
Formal Name: L-leucyl-6-diazo-5-oxo-L-norleucine, ethyl ester
MF: $C_{14}H_{24}N_4O_4$
FW: 312.4
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: -80°C
Stability: ≥ 4 years



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

Laboratory Procedures

JHU083 is supplied as a crystalline solid. A stock solution may be made by dissolving the JHU083 in the solvent of choice, which should be purged with an inert gas. JHU083 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

JHU083 is a prodrug form of the glutaminase inhibitor 6-diazo-5-oxo-L-nor-leucine (Item No. 17580).¹ It decreases tumor volume, increases tumor microenvironment (TME) levels of Tnf-secreting macrophages, and increases survival in a B6CaP murine prostate cancer model.¹ JHU083 increases tumor infiltration by Cd8⁺ T cells in a UN-SCC680 murine lung squamous cell carcinoma model.² It also reduces granuloma area and lung weight and increases survival in *M. tuberculosis*-infected mice.³

References

1. Praharaj, M., Shen, F., Lee, A.J., *et al.* Metabolic reprogramming of tumor-associated macrophages using glutamine antagonist JHU083 drives tumor immunity in myeloid-rich prostate and bladder cancers. *Cancer Immunol. Res.* **12**(7), 854-875 (2024).
2. Huang, M., Xiong, D., Pan, J., *et al.* Targeting glutamine metabolism to enhance immunoprevention of EGFR-driven lung cancer. *Adv. Sci. (Weinh)* **9**(26), e2105885 (2022).
3. Parveen, S., Shen, J., Lun, S., *et al.* Glutamine metabolism inhibition has dual immunomodulatory and antibacterial activities against *Mycobacterium tuberculosis*. *Nat. Commun.* **14**(1), 7427 (2023).

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, 05/11/2026

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