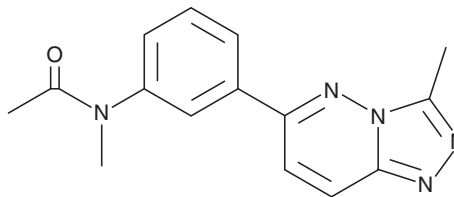


# PRODUCT INFORMATION



**Lin28 1632**  
Item No. 22401

**CAS Registry No.:** 108825-65-6  
**Formal Name:** N-methyl-N-[3-(3-methyl-1,2,4-triazolo[4,3-b]pyridazin-6-yl)phenyl]-acetamide  
**MF:** C<sub>15</sub>H<sub>15</sub>N<sub>5</sub>O  
**FW:** 281.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 247, 285 nm  
**Supplied as:** A crystalline 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

Lin28 1632 is supplied as a crystalline solid. A stock solution may be made by dissolving the lin28 1632 in the solvent of choice, which should be purged with an inert gas. Lin28 1632 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of lin28 1632 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of lin28 1632 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of lin28 1632 in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

## Description

Lin28 1632 is a small molecule inhibitor of the interaction between the RNA binding protein Lin28 and let-7 precursor RNA (IC<sub>50</sub> = 8 μM in a competition ELISA).<sup>1</sup> It increases levels of endogenous let-7 miRNAs and decreases Lin28 expression in a dose-dependent manner in murine embryonic stem cells (mESCs). Lin28 1632 inhibits clonogenic growth of 22Rv1, PC3, DU145, and Huh7 cancer cells and decreases tumor-sphere formation by 22Rv1 and Huh7 cells *in vitro*. It also binds bromodomain-containing protein 4 (BRD4) and CREB-binding protein (CBP/CREBBP) bromodomains (K<sub>d</sub>s = 7 and 25 μM, respectively).

## Reference

1. Roos, M., Pradère, U., Ngondo, R.P., *et al.* A small-molecule inhibitor of Lin28. *ACS Chem Biol.* **11**(10), 2773-2781 (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, 11/28/2022

## 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