

# PRODUCT INFORMATION



## HeLa Whole Cell Lysate

Item No. 31640

### Overview and Properties

---

**Source:** Human cervix epitheloid carcinoma cell line  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Supplied in:** Lysed in diluted RIPA Buffer (Item No. 10010263) including protease inhibitors  
**Concentration:** *batch specific* mg/ml  
*Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.*

### Description

---

HeLa whole cell lysate is intended for use as a control for various biological assays.

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, 10/28/2020

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