

PRODUCT INFORMATION



K562 Whole Cell Lysate

Item No. 31741

Overview and Properties

Source: Human myelogenous erythroleukemia 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

K562 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, 10/23/2023

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