

# Manual for 5x Lysis Buffer



Trade name: **5x Universal Lysis Buffer**

Product #: 333

---

## Product description

Product name: 5x Universal Lysis Buffer

Product No.: 333

Manufacturer: Prolume / NanoLight

Quantity: 50 ml

Buffer systems: Use with the following buffer systems:

Cat. # 318 Firefly Assay Reagent

Cat. # 324 NLuc FLASH buffer

Storage conditions: Long term storage at -20°C. Short term (several months) at 4°C. Let product come to room temperature before use.

## Manual

1. Although this is a 5x concentrated lysis buffer, we recommend adjusting the concentration according to your cell line system. See the table below:

Cell Line	Recommended Lysisbuffer dilution	Notes
HEK293/HEK293T	20x to 5x	10x for cytosolic lysis, 5x for stronger solubilization
HeLa	10x to 5x	5x for full protein lysis
CHO	10x to 5x	Usually using the buffer as 10x works very well
NIH 3T3	5x	Fibroblasts, use as 5x
Jurkat / K562	20x	Very susceptible to lysis use less lysis buffer
C2C12 (muscle)	5x to 2x	Tougher membrane, combine with freeze-thaw
Primary cells	20x to 10x	Start with lower amounts first

2. For a standard lysis solution prepare 1x lysis buffer by adding 4 volumes of double distilled water to 1 volume of 5x lysis reagent. For a milder lysis mix 9 volumes of double distilled water and 1 volume of 5x lysis reagent. Add to washed cells as followed.

# ***Manual for 5x Lysis Buffer***



Trade name: **5x Universal Lysis Buffer**

Product #: 333

---

## **Preparing Mammalian Cell lysates:**

1. Remove growth medium from cultured cells. Usually, high amounts of FCS will cause background luminescence with Coelenterazine based assays.
2. Carefully rinse cells with PBS or TBS. Remove wash solution completely.
3. Add the prepared 1x lysis reagent to your culture vessel to cover its surface. (e.g. 20  $\mu$ l/well in 96 well plate, 400  $\mu$ l per 6 cm culture dish, 900  $\mu$ l for 10 cm culture dish)
4. For culture dishes, scrape adherend cells from dish and transfer into microcentrifuge tube. Pellet debris by brief centrifugation and transfer supernatant into a new tube.
5. Follow the instructions of your luciferase assay kit. In general mix equal amounts of lysate (e.g. 50  $\mu$ l) with the luciferase assay buffer (e.g. 50  $\mu$ l FLuc buffer + added D-Luciferin from the #318 kit).