

ab65393 LDH-Cytotoxicity Assay Kit II

For the rapid, sensitive and accurate measurement of LDH-Cytotoxicity in cell culture samples.
This product is for research use only and is not intended for diagnostic use

Storage and Stability: All components in this kit are shipped on blue ice and are suitable for storage at -20°C, unless reconstituted. Upon receipt, immediately store kit at -20°C in the dark. Individual components may be stored at alternative temperatures as shown in the table below. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Materials Supplied:

| Item | Quantity | Storage Temperature (Before Prep) |
|----------------------|----------|-----------------------------------|
| Substrate Mix | 1 vial | -20°C |
| LDH Assay Buffer | 50 mL | -20°C |
| Lysis Buffer II | 5 mL | -20°C or Ambient |
| Stop Solution IV | 5 mL | -20°C |
| LDH Positive Control | 1 vial | -20°C |

Materials Required, Not Supplied

- These materials are not included in the kit, but will be required to successfully perform this assay:
- Microcentrifuge
- Pipettes and pipette tips
- Colorimetric microplate reader
- 96 well plate
- Orbital shaker

Assay Protocol

1. Reconstitute LDH Positive Control with 100 µl of LDH Assay Buffer.
Note: Aliquot LDH Positive Control after reconstitution. Store aliquots at -20 °C. Avoid freeze/thaw cycles.
2. **Sample Preparation:** Collect cells (adherent or suspension) and wash once with fresh regular culture medium, then seed 100 µl cells (with 2-10 x 10⁴ cells*) in a 96-well plate as the following:
 - a. **Background Control:**
100 µl culture medium per well in triplicates with no cells. The Background Control will measure reagents and LDH background from culture medium serum. The background value has to be subtracted from all other values.
 - b. **Low Control:**
100 µl cells in triplicate wells
 - c. **High Control:**

100 µl cells in triplicates, add 10 µl Lysis Buffer II each well, mix. To adjust the increase of medium volume, 11 µl of the medium may be used in LDH activity assay at step 5.

d. Test Sample:

100 µl cells in triplicates, add test substances each well, mix.

Notes:

- a) Trypsin may be used to remove adherent cells from a culture surface before seeding in a 96-well plate.
 - b) The amount of cells to be used per well depends on the cell types. To optimize the assay, you can do a quick testing by using 2, 4, 8 x 10⁴ cells per well, and then follow the assay protocol to determine the cell number you should use. The high control should be OD_{450nm} ~2.0 after 30 min treatment with 10 % Lysis Buffer II, while the low control should be OD_{450nm} < 0.8. The reaction time should be set at ~ 30 min.
 - c) LDH Positive Control (5 µl LDH) can be used to test whether all reagents are working properly to response to active LDH enzyme.
 - d) If the test substances are not dissolved in PBS, a solvent control may be performed by addition of the same amount of solvent in triplicates without testing substances.
3. **Sample Incubation:** Incubate cells in an incubator (5 % CO₂, 90 % humidity, 37°C) for the appropriate time of treatment determined for test substance. Gently shake the plate at end of the incubation to ensure LDH is evenly distributed in the culture medium.
 4. Centrifuge cells at 600 x g for 10 min to precipitate the cells.
 5. Transfer the clear medium solution (10 µl/well) into an optically clear 96-well plate.
 6. **LDH Reaction Mix:** Reconstitute the Substrate Mix in 1.1 ml ddH₂O for 10 min and mix thoroughly. The solution is stable for two months at 4°C.
For 100 assays, mix 200 µl of Substrate Mix with 10.0 ml of LDH Assay Buffer. The LDH Reaction Mix should be stable for several weeks at 4°C. Add 100 µl LDH Reaction Mix to each well, mix and incubate for 30 min** at room temperature.
 7. Measure the absorbance of all controls and samples with a plate reader equipped with 450 nm (440 - 490 nm) filter. The reference wavelength should be 650 nm.

Notes:

- a) The reaction time can be decreased or increased depending on the color development. The plate can be read at multiple time points until the desired reading is observed. The high control should be OD_{450nm} ~2.0, while the low control should be OD_{450nm} < 0.8.
- b) The reaction can be stopped by adding 10 µl of Stop Solution IV, mix and read within 48 hours without significant changes. Protect the reaction from light and evaporation.

Data analysis

$$\text{Cytotoxicity (\%)} = \frac{\text{(Test Sample - Low Control)}}{\text{(High Control - Low Control)}} \times 100$$

Troubleshooting

| Problem | Reason | Solution |
|------------------------------------|---|--|
| Assay not working | Assay buffer at wrong temperature | Assay buffer must not be chilled - needs to be at RT |
| | Protocol step missed | Re-read and follow the protocol exactly |
| | Plate read at incorrect wavelength | Ensure you are using appropriate reader and filter settings (refer to datasheet) |
| | Unsuitable microtiter plate for assay | Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells |
| Unexpected results | Measured at wrong wavelength | Use appropriate reader and filter settings described in datasheet |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Unsuitable sample type | Use recommended samples types as listed on the datasheet |
| | Sample readings are outside linear range | Concentrate/ dilute samples to be in linear range |
| Samples with inconsistent readings | Unsuitable sample type | Refer to datasheet for details about incompatible samples |
| | Samples prepared in the wrong buffer | Use the assay buffer provided (or refer to datasheet for instructions) |
| | Samples not deproteinized (if indicated on datasheet) | Use the 10kDa spin column (ab93349) |
| | Cell/ tissue samples not sufficiently homogenized | Increase sonication time/ number of strokes with the Dounce homogenizer |
| | Too many freeze-thaw cycles | Aliquot samples to reduce the number of freeze-thaw cycles |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Samples are too old or incorrectly stored | Use freshly made samples and store at recommended temperature until use |

| Problem | Reason | Solution |
|---|--|--|
| Standard curve is not linear | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| | Pipetting errors when setting up the standard curve | Try not to pipette too small volumes |
| | Incorrect pipetting when preparing the reaction mix | Always prepare a master mix |
| | Air bubbles in wells | Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates |
| | Concentration of standard stock incorrect | Recheck datasheet for recommended concentrations of standard stocks |
| | Errors in standard curve calculations | Refer to datasheet and re-check the calculations |
| Lower/ Higher readings in samples and standards | Use of other reagents than those provided with the kit | Use fresh components from the same kit |
| | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| | Out-of-date kit or incorrectly stored reagents | Always check expiry date and store kit components as recommended on the datasheet |
| | Reagents sitting for extended periods on ice | Try to prepare a fresh reaction mix prior to each use |
| | Incorrect incubation time/ temperature | Refer to datasheet for recommended incubation time and/ or temperature |
| | Incorrect amounts used | Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume) |

Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)