

Cytotoxicity LDH Assay Kit-WST

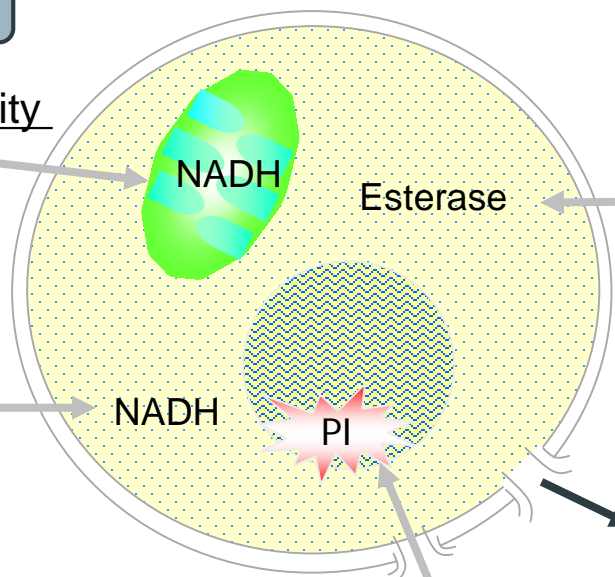
Cell proliferation assay / Cell death assay

Cell proliferation assay

Mitochondrial respiration activity

Dehydrogenase activity

Cell Counting Kit -8



Esterase activity

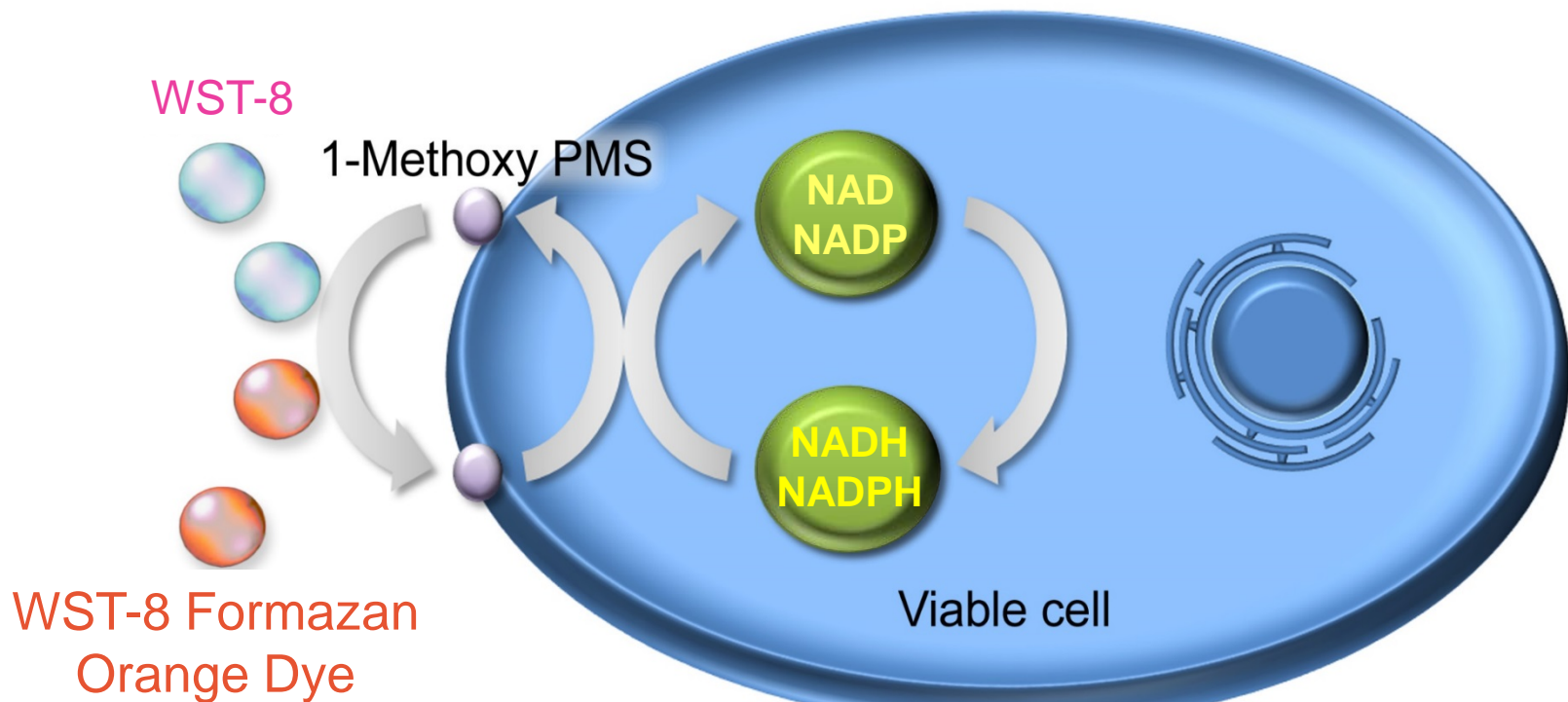
Cell death assay

Released LDH activity

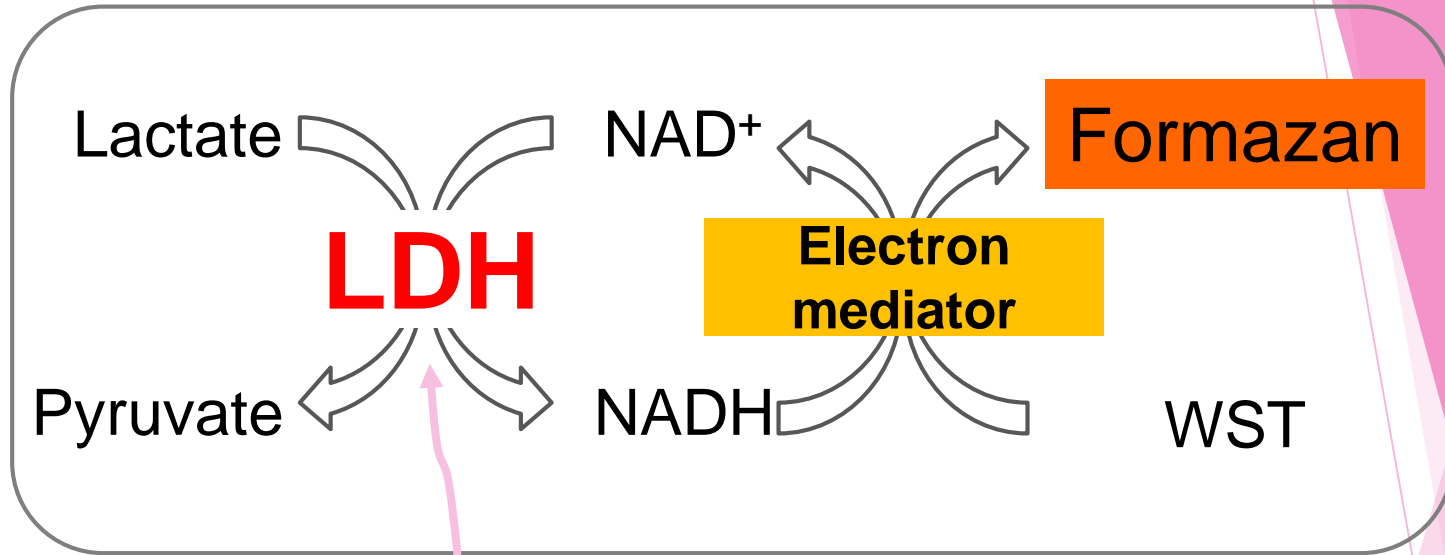
Cytotoxicity LDH Assay Kit-WST

Cell membrane damage

Principle of Cell Counting Kit-8 Assay



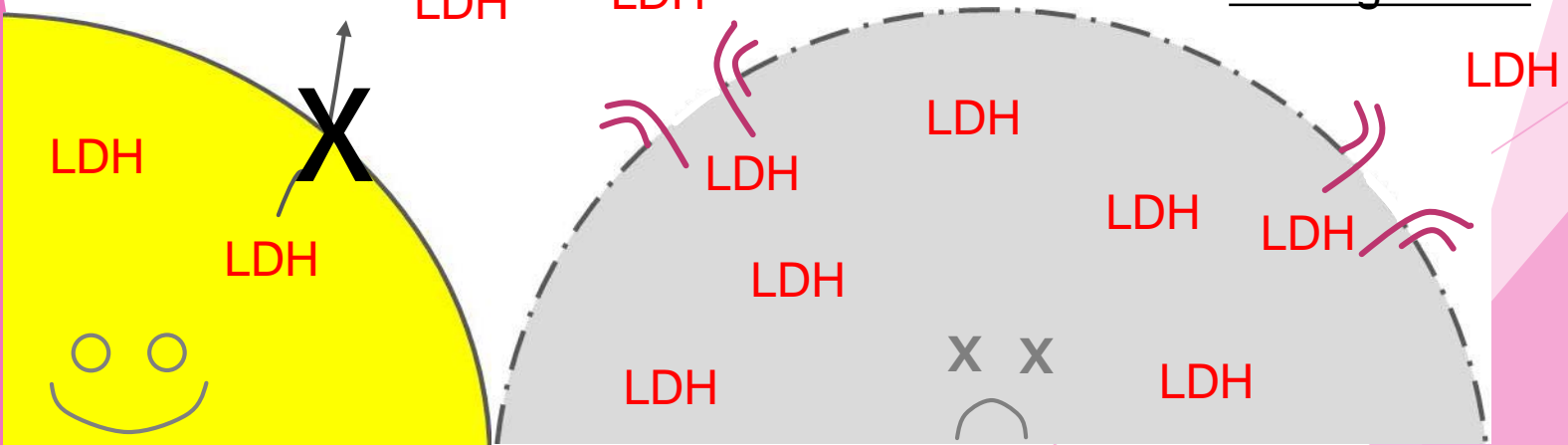
Principle of LDH-release cytotoxicity assay



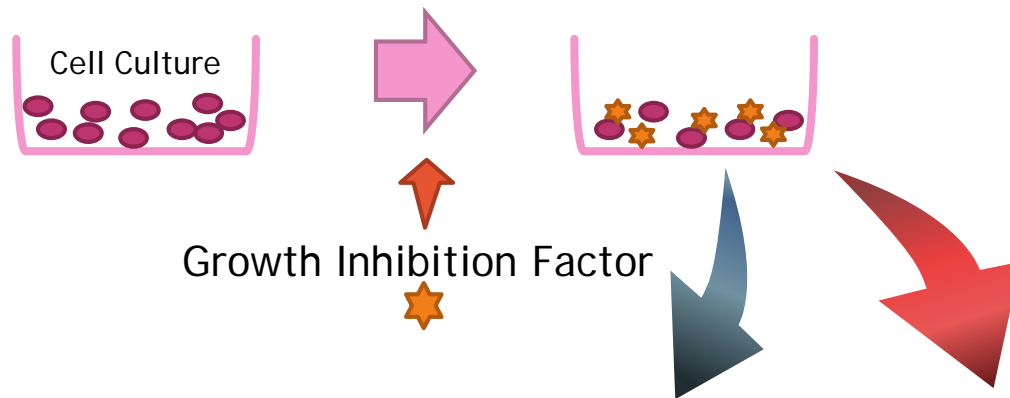
Live Cell

LDH-release

Damaged Cell

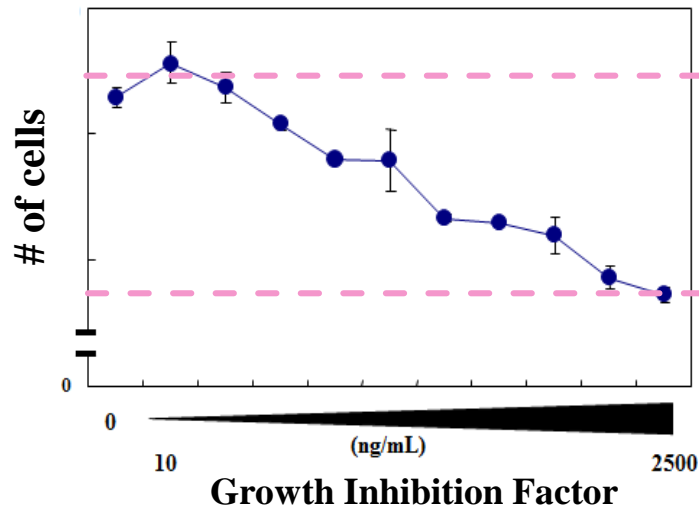


Application (For Example)



Growth Inhibition Factor

Cell Counting Kit-8 Assay



Cytotoxicity LDH Assay Kit-WST

*Effect from the factor?
or
Damage to the cell?*

Cytotoxicity LDH Assay Kit-WST for Cell Death Assay

Feature

- ✓ Both Homogeneous and Non-homogeneous Assays Possible
- ✓ Longer Working Solution Stability(0-5°C for 2 months)
No need to freeze the working solution
- ✓ High Linearity

Kit Contents

- Dye Mixture ×1
- Assay Buffer ×1
- Lysis Buffer ×1
- Stop Solution ×1

Storage condition

- Stored at 0 - 5 °C

Stability

- Stable at 0 - 5 °C for one year



Cytotoxicity LDH Assay Kit-WST

Both Homogeneous and Non-homogeneous Assay Possible

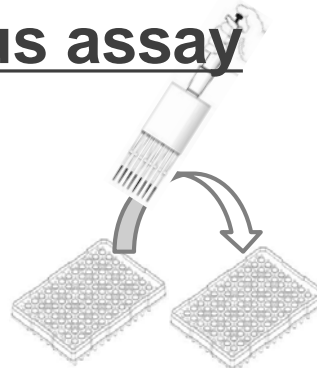
Homogeneous assay



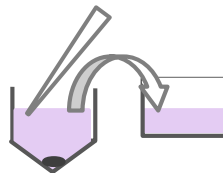
Working Solution



Non-homogeneous assay



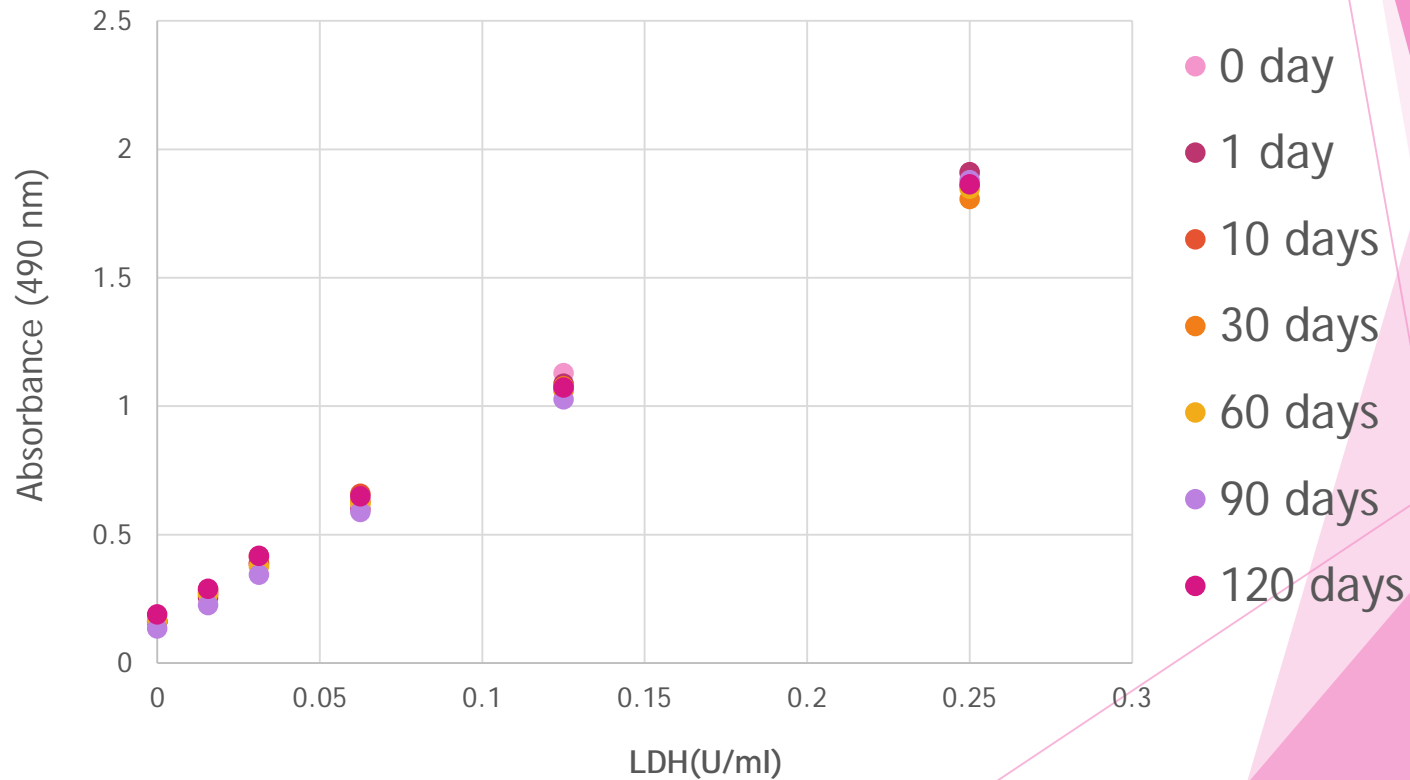
Working Solution



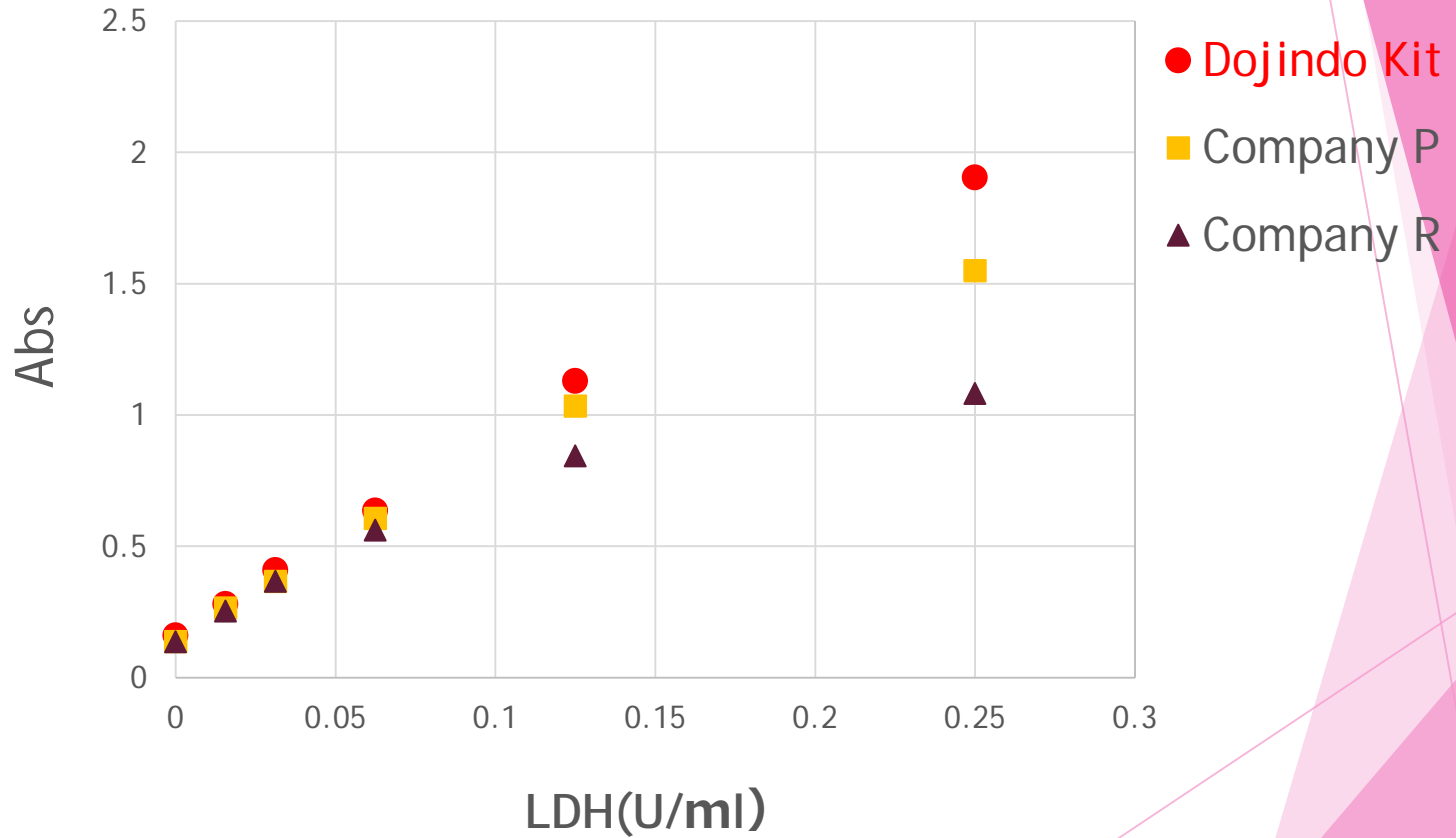
Longer Working Solution Stability

Storage Condition: 0-5 °C for 2 months

Cytotoxicity LDH Assay Kit-WST



High Linearity



Price

| | |
|-------------|------------------------|
| 100 tests: | \$80.00 |
| 500 tests: | \$198.00 |
| 2000 tests: | \$398.00 (Coming soon) |

Cytotoxicity LDH Assay Kit-WST for Cell Death Assay

Feature

- ✓ Both homogeneous and non-homogeneous assays possible
- ✓ Longer Working solution stability(0-5°C for 2 months)
No need to freeze the working solution
- ✓ High Linearity

✓ Detail Information of
"Cytotoxicity LDH Assay Kit-WST"

✓ Promo Here!!



Cytotoxicity LDH Assay Kit-WST