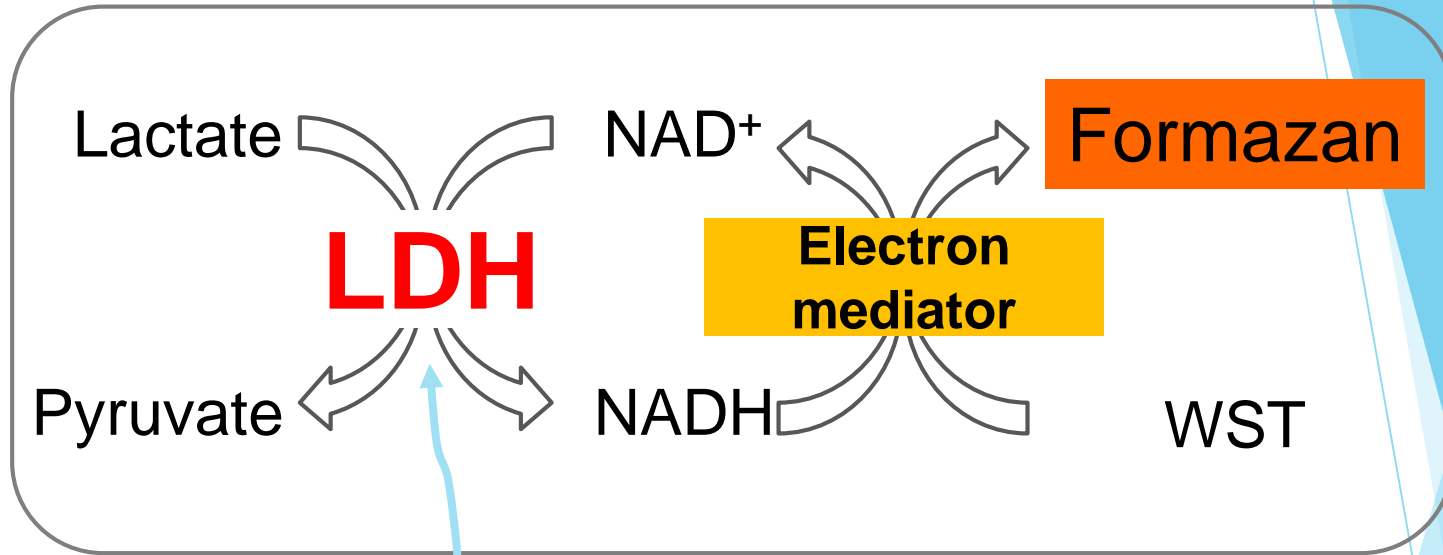


Cytotoxicity LDH Assay Kit-WST

Principle of LDH-release cytotoxicity assay



Live Cell

LDH-release

Damaged Cell

LDH

LDH

LDH

LDH

LDH

LDH

LDH

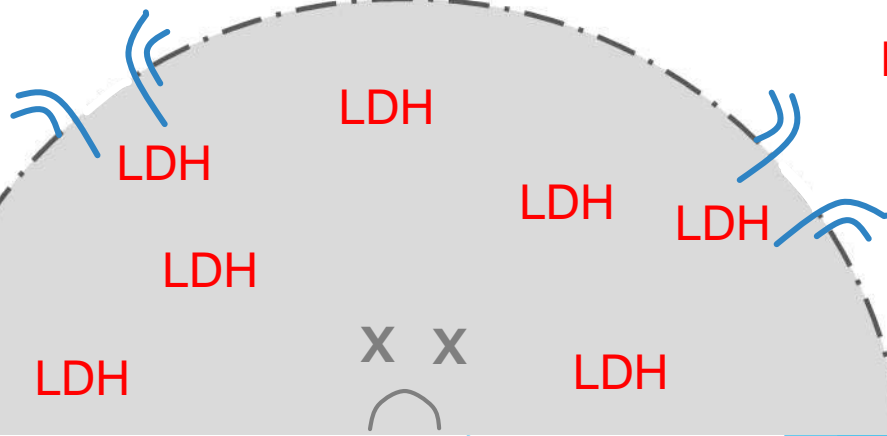
LDH

LDH

LDH

LDH

LDH



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

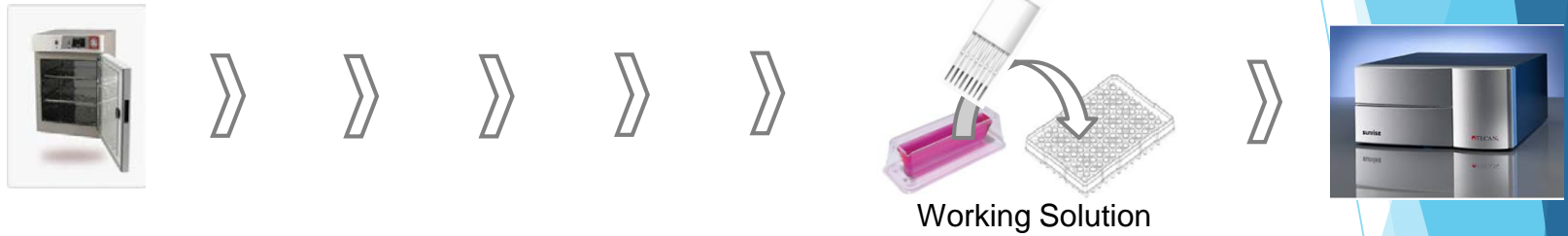
- Stable at 0 - 5 °C for one year



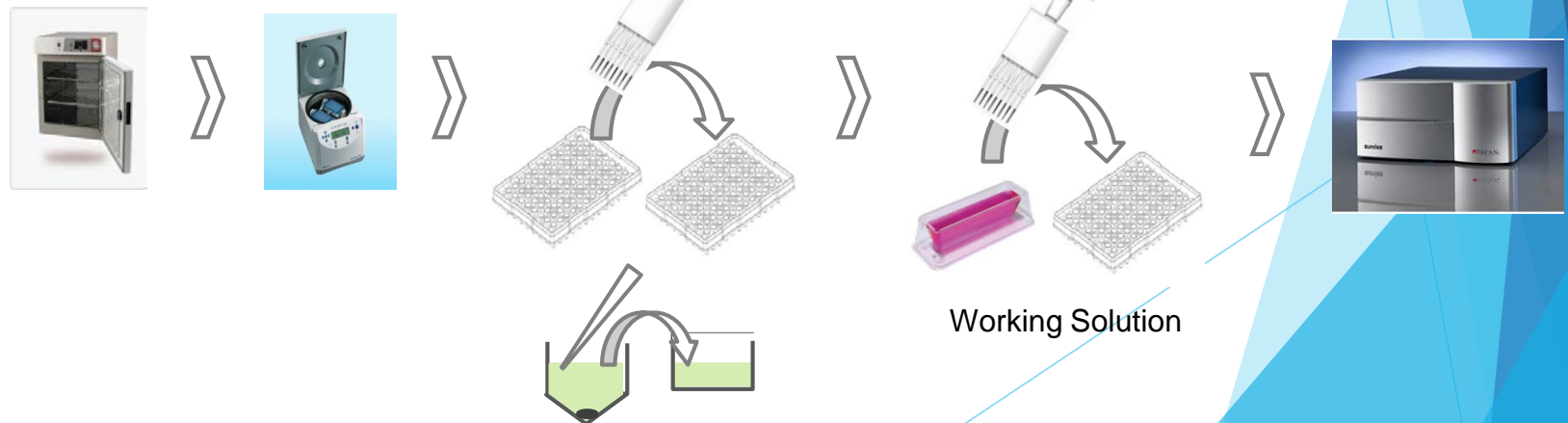
Cytotoxicity LDH Assay Kit-WST

Both Homogeneous and Non-homogeneous Assay Possible

Homogeneous assay



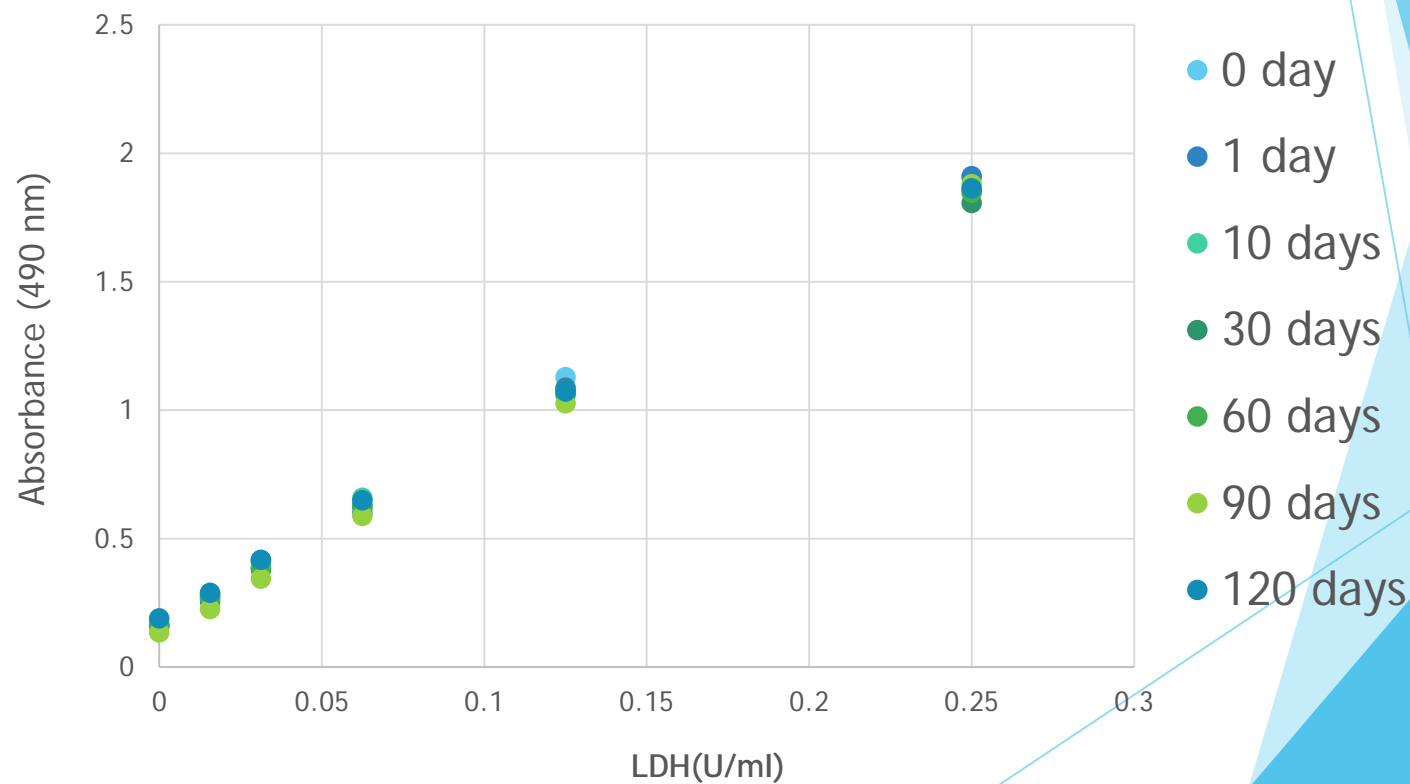
Non-homogeneous assay



Longer Working Solution Stability

Storage Condition: 0-5 °C for 2 months

Cytotoxicity LDH Assay Kit-WST



Longer Working Solution Stability

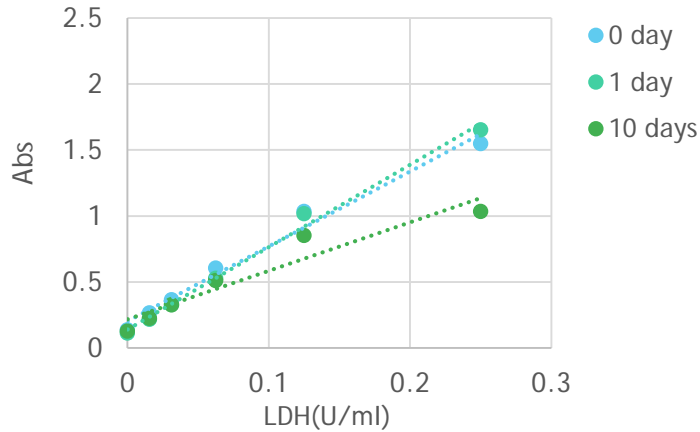
Product

Company A
Company B
Company B`
Company C

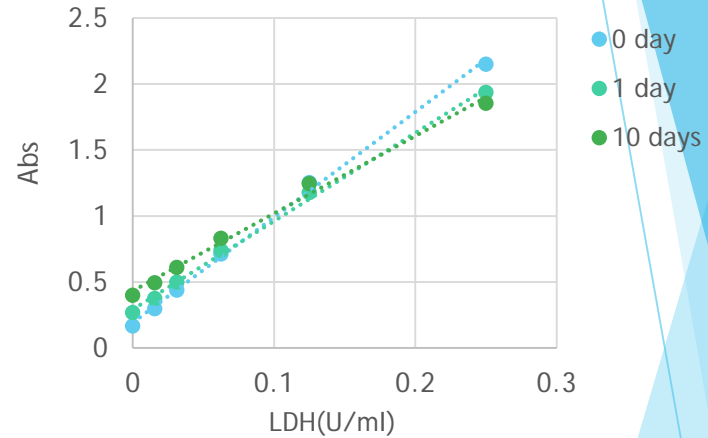
Strage Condition

-20 °C for 6-8 weeks
5 °C for several weeks
×(N/A)
×(N/A)

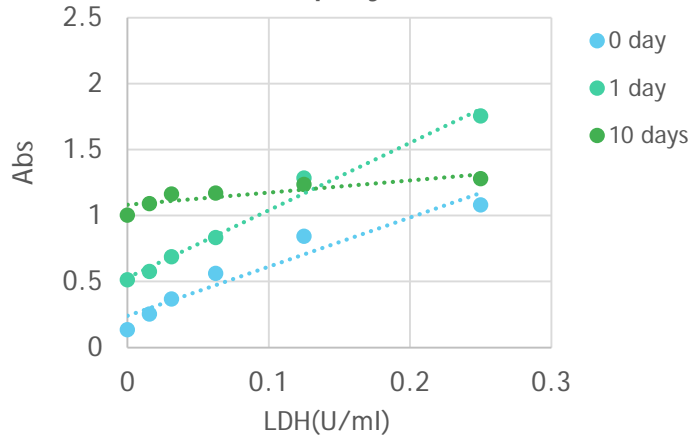
Company A



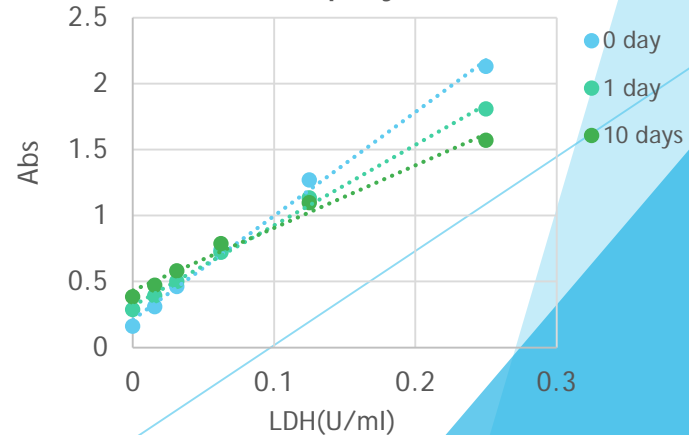
Company B



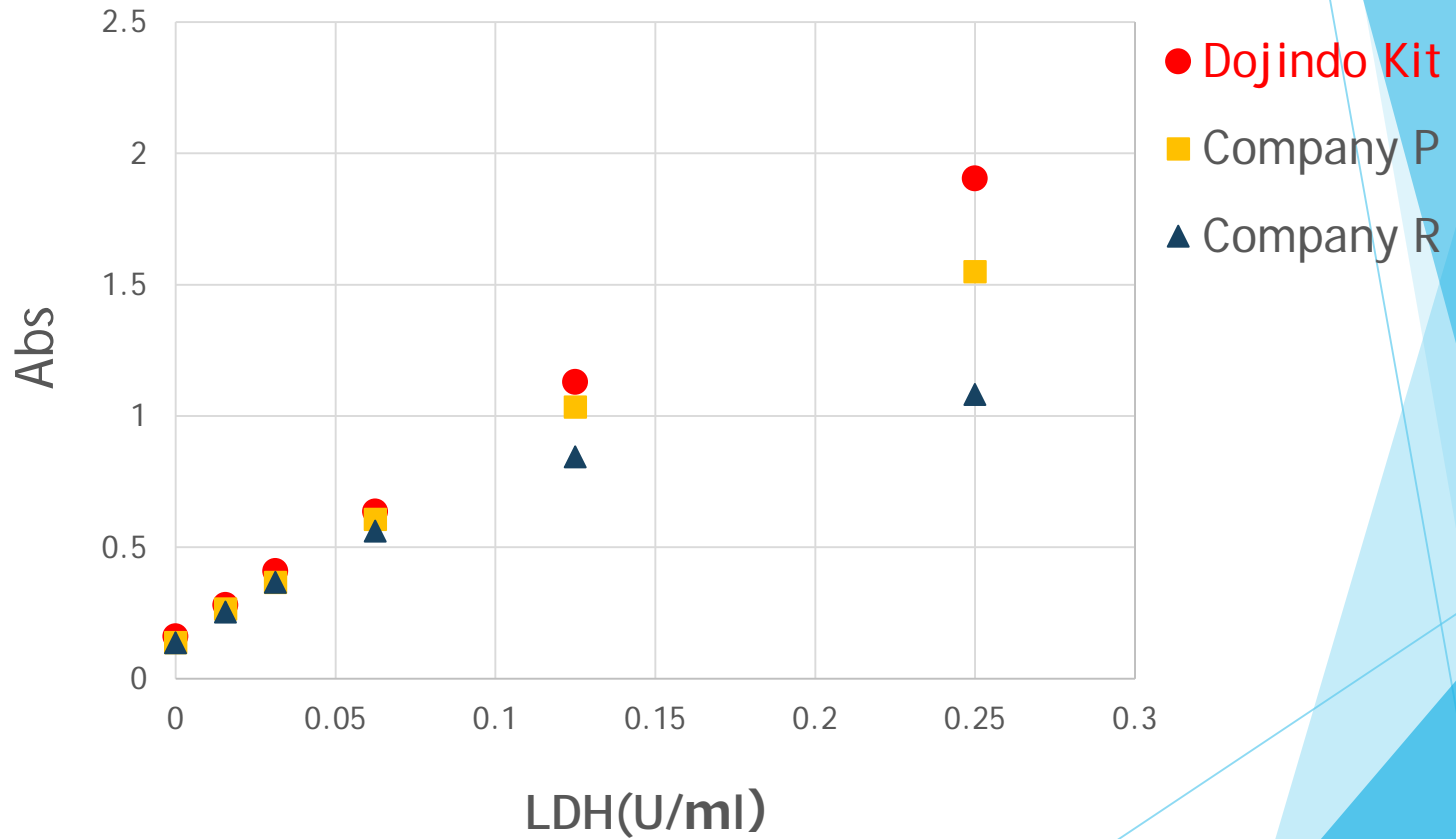
Company B`



Company C



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