

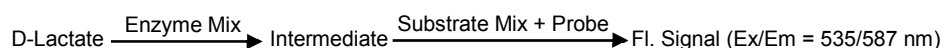
PicoProbe™ D-Lactate Fluorometric Assay Kit

rev. 4/13

(Catalog # K668-100; 100 assays; Store at -20°C)

I. Introduction:

D-Lactate is the result of anaerobic glycolysis by microorganisms in the gastrointestinal system, and the product of detoxification of methylglyoxal by the glyoxalase system. The presence of abnormal levels of D-Lactate has been linked to a series of pathological conditions, such as diabetes, and appendicitis. BioVision's PicoProbe™ D-Lactate Assay kit offers simplicity, sensitivity, and can be adapted to high-throughput research. The assay enzymatically oxidizes D-Lactate generating a fluorescent signal (Ex/Em = 535/587 nm). The signal is directly proportional to the amount of D-Lactate. The assay kit can detect D-Lactate as low as 0.1 μM in a variety of samples.



II. Application:

- Measurement of D-Lactate in various tissues/cells/biological fluids
- Analysis of D-lactate in pathological conditions
- Mechanistic study of the glyoxalase system

III. Sample Type:

- Serum, plasma, urine & other body fluids
- Animal tissues: liver, muscle, heart, etc.
- Cell culture: adherent or suspension cells
- Fermentation media
- Food

IV. Kit Contents:

Components	K668-100	Cap Code	Part Number
D-Lactate Assay Buffer	25 ml	WM	K668-100-1
PicoProbe™ (in DMSO)	0.4 ml	Blue	K668-100-2
D-Lactate Enzyme Mix (Lyophilized)	1 vial	Green	K668-100-3
D-Lactate Substrate Mix (Lyophilized)	1 vial	Red	K668-100-4
D-Lactate Standard (100 mM)	0.1 ml	Yellow	K668-100-5

V. User Supplied Reagents and Equipment:

- White 96-well plate with flat bottom
- Multi-well spectrophotometer

VI. Storage and Handling:

Store kit at -20°C, protected from light. Warm D-Lactate Assay Buffer to room temperature before use. Briefly centrifuge small vials prior to opening.

VII. Reagent Preparation and Storage Conditions:

- **PicoProbe™:** Ready to use as supplied. Warm to room temperature before use. Store at -20°C.
- **D-Lactate Enzyme Mix:** Reconstitute with 220 μl D-Lactate Assay Buffer. Pipette up and down to dissolve completely. Aliquot and store at -20°C. Avoid repeated freeze/thaw. Keep on ice while in use. Stable for 2 months at -20°C.
- **D-Lactate Substrate Mix:** Reconstitute with 220 μl D-Lactate Assay Buffer. Pipette up and down to dissolve completely. Aliquot and store at -20°C. Avoid repeated freeze/thaw. Keep on ice while in use. Stable for 2 months at -20°C.

VIII. D-Lactate Assay Protocol:

1. **Sample Preparation:** Prepare 2-50 μl test samples in a 96-well plate (for high target concentrations, dilute the samples). Bring the volume to 50 μl/well with D-Lactate Assay Buffer. We suggest using different volumes of your sample to ensure the readings are within the Standard Curve range.

Notes:

- Tissue (20 mg) or cells (2×10^6) can be homogenized in 100 μl D-Lactate Assay Buffer. Centrifuge at 10,000x g for 10 min. to remove insoluble material. Soluble fractions may be assayed directly.
 - Body fluids, such as urine, should be filtered using a 10 kDa MW spin filter (BioVision, Cat # 1997-25). Dilute samples with pure water. An appropriate dilution factor is usually between 1:10 and 1:100. Diluted samples can be assayed directly.
 - Food samples. A) Beer: Remove CO₂ by vacuum-filtering samples. Dilute samples with pure water (appropriate dilution factor is approximately 1:10). Diluted samples can be assayed directly. B) Yogurt: vortex 1 gram of yogurt in 10 ml of water until homogeneous. Dilute with pure water (dilution factor ~ 1:100 to 1:500). Diluted samples can be assayed directly.
 - Endogenous enzyme activity may cause loss of D-Lactate. Samples containing enzyme activity (such as culture medium or tissue lysate) should be kept at -80°C or filtered through a 10 kDa MW spin filter (BioVision, Cat.# 1997-25).
2. **Standard Curve Preparation:** Dilute D-Lactate Standard to 1 mM (1000 pmol/μl) by adding 10 μl of 100 mM D-Lactate Standard to 990 μl D-Lactate Assay Buffer, mix well. Dilute further to 0.02 mM (20 pmol/μl) by adding 20 μl of 1 mM D-Lactate Standard to 980 μl of D-

Lactate Assay Buffer. Mix well. Add 0, 2, 4, 6, 8 & 10 μl of 0.02 mM D-Lactate Standard into a series of wells in a 96 well plate to generate 0, 40, 80, 120, 160 and 200 pmol/well of D-Lactate Standard. Adjust volume to 50 μl /well with D-Lactate Assay Buffer.

- 3. Reaction Mix:** Mix enough reagents for the number of assays (samples and Standards) to be performed. For each well, prepare 50 μl Reaction Mix containing:

	Reaction Mix	Background Control Mix*
D-Lactate Assay Buffer	45 μl	47 μl
PicoProbe™	1 μl	1 μl
D-Lactate Enzyme Mix	2 μl	---
D-Lactate Substrate Mix	2 μl	2 μl

Add 50 μl of the Reaction Mix to each well containing Standard & test samples. Mix well.

***Note:** For samples having high NADH levels, add 50 μl of Background Control Mix to sample background control well(s). Mix well.

- 4. Measurement:** Incubate the reaction for 30 minutes at 37°C, protected from light. Measure fluorescence (Ex/Em = 535/587 nm) in a microplate reader with kinetics mode.
- 5. Calculation:** Subtract 0 D-Lactate Standard reading from all readings. Plot the D-Lactate Standard curve. If sample background control reading is significantly high, subtract the background control reading from sample reading. Apply the corrected sample reading to the D-Lactate Standard curve to get B pmol of D-Lactate in the sample wells.

$$\text{Sample D-Lactate concentration} = B/V \times \text{Dilution Factor} = \text{pmol}/\mu\text{l} = \text{nmol}/\text{ml} \text{ or } \mu\text{M}$$

Where: **B** = amount of D-Lactate in sample well from Standard curve (pmol)

V = sample volume added in the reaction well (μl)

D-Lactate in samples can also be expressed in nmol/mg of sample.

D-Lactate molecular weight: 112.1 g/mol.

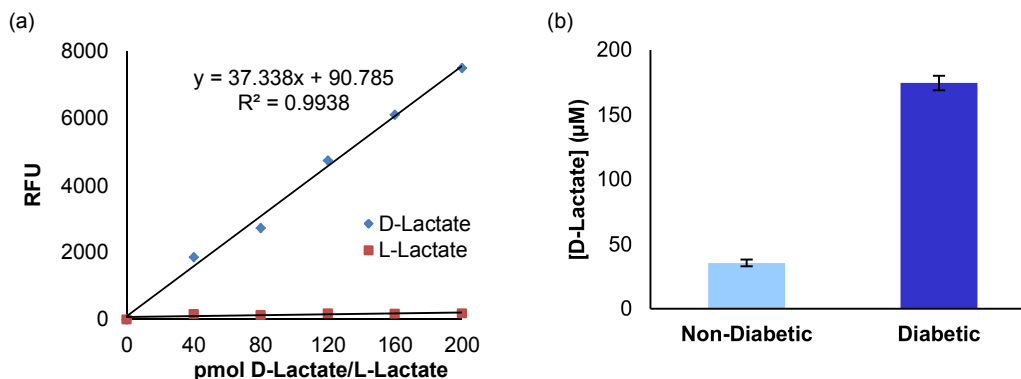


Figure: (a) D-Lactate Standard Curve. The assay specifically detects D-Lactate, not L-Lactate. (b) Measurement of D-Lactate in urine: Diluted samples (1:10, non-diabetic; 1:100 diabetic) were spiked with known amounts of D-Lactate (0-200 pmol) and assayed as specified.

IX. RELATED PRODUCTS:

Lactate Colorimetric Assay Kit II
 PicoProbe™ Lactate Fluorometric Assay Kit
 Lactate Dehydrogenase (LDH) Activity Assay Kit
 Human Recombinant LDHA
 LDH Antibody
 Glutathione Assay Kit
 Cytotoxicity Assay Kit
 Creatinine Assay Kit
 NADH/NADPH Assay Kits

Lactate Colorimetric/Fluorometric Assay Kit
 D-Lactate Colorimetric Assay Kit
 LDH-Cytotoxicity Colorimetric Assay Kit II
 Human Recombinant LDHB
 Cholesterol Assay Kit
 Glucose Assay Kit II
 Pyruvate Assay Kit
 Cell Proliferation Assays

FOR RESEARCH USE ONLY! Not to be used on humans.