

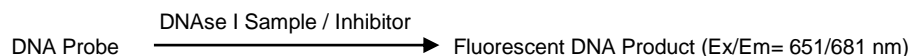
## DNase I Activity Assay Kit (Fluorometric)

3/17

(Catalog # K429-100; 100 assays; Store at -20°C)

### I. Introduction:

Deoxyribonuclease I (DNase I) is an endonuclease that cleaves DNA phosphodiester bonds yielding 5'-phosphorylated and 3'-hydroxylated oligonucleotides. DNase I targets single-stranded DNA, double-stranded DNA, and chromatin in a non-specific manner. As an important player in cellular waste management, DNase I is normally secreted extracellularly to clear the system from circulating cell-free DNA, foreign DNA from food digestion or potential pathogens, and endogenous chromosomal DNA from apoptotic and necrotic cells. Abnormal DNase I activity occurs in association with a variety of cancers and auto-immune illnesses that exhibit elevated levels of cell-free DNA. Furthermore, DNase I has been therapeutically used in cystic fibrosis patients to degrade DNA and reduce sputum viscosity. BioVision's DNase I Activity Fluorometric Assay Kit allows for quantitative evaluation of DNase I activity of purified enzymes and their inhibitors as well as comparative examination of DNase I activity in biological samples. Enzyme activity is detected upon cleavage of a DNA Probe, which yields a fluorescent DNA product measured at Ex/Em = 651/681 nm. The limit of quantification (L.O.Q) is 178 fmoles of DNA probe cleaved per minute per ml.



### II. Applications:

- Measurement of DNase I activity of purified proteins
- Quantitative analysis of DNase I mutants and inhibitors
- Comparative examination of DNase I activity in biological samples

### III. Sample Type:

- Purified Protein, serum.

### IV. Kit Contents:

Components	K429-100	Cap Code	Part Number
10X DNase I Assay Buffer	1.1 ml	Clear	K429-100-1
DNA Probe	1 vial	Blue	K429-100-2
DNA Probe Re-suspension Buffer	250 µl	Red	K429-100-3
DNase I Positive Control	1 vial	Purple	K429-100-4
Positive Control Re-suspension Buffer	1 ml	Brown	K429-100-5
Molecular Biology Grade Water	25 ml	NM	K429-100-6

### V. User Supplied Reagents and Equipment:

- 96-well white plate with flat bottom, low-medium binding
- Spectrophotometer
- Purified DNase I, DNase I inhibitors, biological samples

### VI. Storage Conditions and Reagent Preparation:

Store kit at -20°C, protected from light. Briefly centrifuge small vials prior to opening. Upon re-suspension, aliquot DNA Probe and DNase I Positive control to avoid repeated freeze-thaw cycles. Read entire protocol before performing the assay.

- **10X DNase I Assay Buffer:** Store at 4 °C. Warm to 37 °C temperature before use.
- **DNA Probe:** Reconstitute with 220 µl of DNA Probe Re-suspension Buffer. Aliquot and store at -20°C. Avoid multiple freeze-thaw cycles.
- **DNA Probe Re-suspension Buffer:** Ready to use. Store at RT.
- **DNase I Positive Control:** Reconstitute with 220 µl of Positive Control Re-suspension Buffer. Aliquot and store at -20°C.
- **Positive Control Re-suspension Buffer:** Ready to use. Store at -20°C.
- **Molecular Biology Grade Water:** Ready to use. Store at RT.

### VII. DNase I Activity Assay Protocol:

**Caution!** It is imperative to use molecular biology grade water for sample preparation and filter tips for sample pipetting at all times to avoid DNase contamination.

**1. Sample Preparation:** Thaw any purified enzymes and biological samples along with all the provided assay components on ice, unless otherwise stated. Dilute enzymes, inhibitors, and biological samples to a desired concentration with water or their corresponding storage buffer. Add a desired amount of enzyme, inhibitor, or biological sample to each well and adjust the volume to 50 µl with water. Use water only (no enzyme/sample) for background control reaction. For positive control reaction, add 2 µl of DNase I Positive Control to 48 µl of water. Mix well.

#### Notes:

- Do not store enzyme/inhibitor/sample dilutions; discard the dilutions instead.
- The recommended amount of serum sample to use in the assay is 10-25 µl.
- For uncharacterized enzymes, we suggest testing several doses to ensure the reading is within the Standard Curve range.

- d. If the user suspects any non-specific sample DNase activity, 50 mM 2-Nitro-5-thiocyanatobenzoic acid (Cat. # B1316) can be used to specifically inhibit DNase I activity.
2. **DNA Probe Standard Curve:** Prepare 1  $\mu\text{M}$  DNA Probe stock by diluting 4  $\mu\text{l}$  of 25  $\mu\text{M}$  DNA Probe in 96  $\mu\text{l}$  of molecular biology grade water. Add 0, 4, 8, 12, 16, 20  $\mu\text{l}$  of 1  $\mu\text{M}$  DNA Probe into a series of wells on a 96-well plate to generate 0, 4, 8, 12, 16, 20, pmol/well of DNA Probe Standard. Adjust the volume to 50  $\mu\text{l}$  with molecular biology grade water.
3. **Reaction Mix:** Mix enough reagents for the number of assays to be performed. For each well containing sample(s), and standards prepare 50  $\mu\text{l}$  Mix containing:

	<u>Sample Reaction Mix</u> (1 assay)	<u>DNA Probe Standard Reaction Mix</u> (1 assay)
10X DNase I Assay Buffer	10 $\mu\text{l}$	10 $\mu\text{l}$
DNA Probe (25 $\mu\text{M}$ )	2 $\mu\text{l}$	---
DNase I Positive Control	---	2 $\mu\text{l}$
Molecular Biology Grade $\text{H}_2\text{O}$	38 $\mu\text{l}$	38 $\mu\text{l}$

Mix and add 50  $\mu\text{l}$  of the Sample Reaction Mix to each well containing the Positive Control, Test Samples, and Background Control. Add 50  $\mu\text{l}$  of DNA Probe Standard Reaction Mix to each well containing DNA Probe Standard.

4. **Measurement:** For positive control, test samples, background control, and DNA Probe Standard measure fluorescence (Ex/Em = 651/681 nm) in kinetic mode every 30 seconds for at least 90 minutes at 37  $^{\circ}\text{C}$ . Adjust GAIN/PMT setting of your fluorometer as necessary so that the standard curve readings are within the detection range of the instrument.
5. **Calculations:** *Standard Curve:* Record RFU at  $t = 90$  min for each DNA Probe standard curve reading. Plot the DNA Probe standard curve with pmol of DNA on the x-axis and RFU on the y-axis. Apply a linear fit to the DNA standard values and determine the standard curve equation. *Samples/Positive Control:* Apply RFU values at each time point to the standard curve equation to determine pmol of DNA cleaved at each reaction time point. Plot pmol DNA on the y-axis vs. time (in minutes) on the x-axis and determine the slope (pmol/min) of the linear portion of the reaction curve. Subtract background control readings from samples.

$$\text{Sample DNase I Activity} = (\text{slope}/V) \times D \text{ (pmol/min/ml} \equiv \mu\text{U/ml)}$$

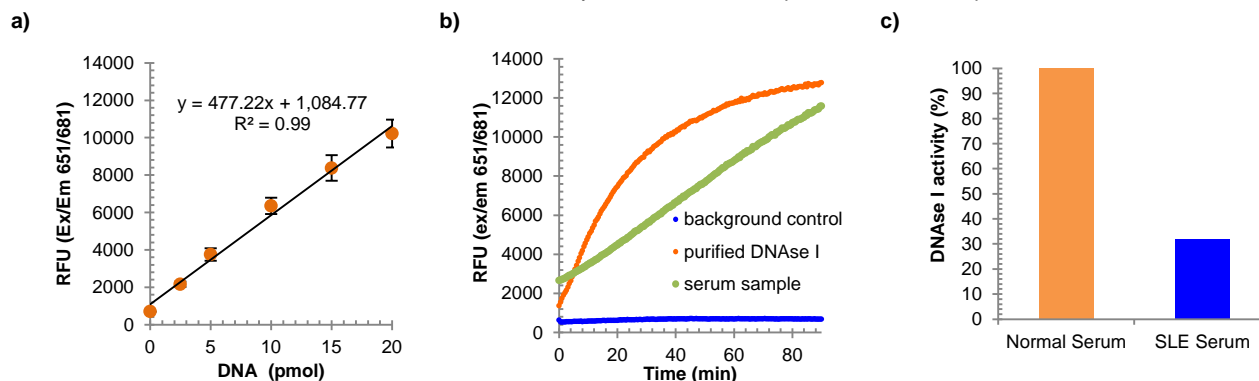
$$\text{Sample Specific Activity} = (\text{slope}/\mu\text{g}) \times D \text{ (pmol/min}/\mu\text{g} \equiv \mu\text{U}/\mu\text{g})$$

Where:  $V$  = sample volume added into the reaction well (ml).

$D$  = Dilution Factor

**Slope** = pmol/min (from the linear range of the activity curve)

**Unit Definition:** One unit of DNase I is the amount of enzyme that cleaved 1.0  $\mu\text{mol}$  of DNA Probe per min. at 37 $^{\circ}\text{C}$ .



**Figure:** a) DNA Probe to Product conversion standard curve; b) representative activity curve for purified DNase I (orange), serum sample (green), and background control (blue) at 37 $^{\circ}\text{C}$ ; c) comparative analysis of DNase I activity from 25  $\mu\text{l}$  undiluted single donor normal vs. Systematic Lupus Erythematosus (SLE) patient serum sample.

## VIII. RELATED PRODUCTS

Apoptotic DNA Ladder Isolation Kit (K170)  
 Quick Apoptotic DNA Ladder Detection Kit (K120)  
 Enhanced Apoptotic DNA Ladder Detection Kit (K130)  
 2-Nitro-5-thiothiocyanatobenzoic acid (B1316)  
 DNase I Antibody (3214)

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