

## Fluoraldehyde™ OPA Protein Assay

### Introduction

The Fluoraldehyde™ o-Phthalaldehyde,(OPA), reagent is a ready-to-use formulation of OPA that offers a rapid fluorescent detection of minute amounts of proteins and peptides in solution. Used in conjunction with the micro-volume capability of the Thermo Scientific NanoDrop™ 3300 Fluorospectrometer, the Fluoraldehyde™ reagent provides a highly sensitive means of protein quantitation with minimal consumption of sample. The ability of the NanoDrop 3300 to measure as little as 2 ul of protein samples allows significantly scaled-down reaction volumes, thereby using only a fraction of sample and reagent commonly needed for conventional cuvette-based fluorometers. **Notes:** there is considerable protein-to-protein and peptide-to-peptide variation with this assay so use of a purified sample of the particular protein or peptide as a standard is recommended. Buffers containing primary amines are not recommended. Peptides or proteins with blocked primary amines will not give a response with OPA.

### Dynamic Range

A wide dynamic range for the assay can be obtained by utilizing two distinct sample to reagent volumes. The high 15 ug/ml to 1000ug/ml range for BSA is covered by mixing 2 ul of protein sample with 20 ul of reagent. The sensitivity of the assay may be increased to 1-50 ug/ml by using a higher sample volume to reagent volume.

### High Range (15 ug/ml–1000 ug/ml)

A 1 to 10 sample to reagent volume is typically used for a high range assay.

### Low range (1 ug/ml– 50 ug/ml)

A 1 to 0.5 sample to reagent volume is typically used for a low range assay.

### Supplies

Equipment:

- NanoDrop 3300 Fluorospectrometer
- 2ul pipettor (low retention tips)

Materials:

- Low lint laboratory wipes
- Amber or foil covered 1.5ml polypropylene tubes or 0.2 ml strip tubes and caps

Reagents:

- Fluoraldehyde™ Reagent, Pierce Cat. # 26025
- Pierce BSA standard cat # 23209 (or other protein/peptide standard)

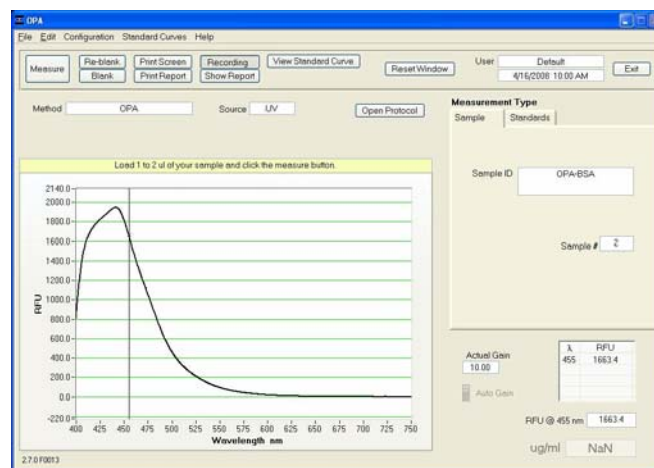
### Assay Recommendations

- Measure 2 ul sample aliquots.
- Mix all solutions gently to avoid micro bubbles.
- Remove samples from the optical surfaces by blotting rather than wiping to reduce residual lint fibers from collecting on the NanoDrop 3300 sampling pedestals.
- Equilibrate all reagents, unknowns and protein standards to room temperature.
- Protect the reagent and solutions containing reagent from light.
- For additional information regarding the Fluoraldehyde™ reagent, please refer to the Pierce web site.

### Fluoraldehyde™ OPA Protein Assay

1. Mix each stock standard solution and unknown samples thoroughly and transfer 10 ul (low range) or 2 ul (high range) of each to the respective amber tubes.
2. Add 5 ul (low range) or 20 ul (high range) of the Fluoraldehyde™ reagent into each standard and sample and mix each standard and unknown sample thoroughly by pipetting up and down several times. Optional: collect the solution at the bottom of the tube by a brief centrifugation,
3. Allow the reaction to incubate 10 minutes at room temperature. Note: It is recommended that the samples be measured immediately after the standard curve has been established.

Example spectrum of Fluoraldehyde™ OPA Protein sample



Rev 4/08

## Standard Curve Protocol

- Clean both sampling pedestals with 2 uL of nuclease free de-ionized water.
- Open upper arm and firmly blot the two pedestals with a dry lab wipe. Make sure there are no traces of lint on the pedestals before continuing.
- Open the operating software. Click on the Protein Quantitation button and select the Flouraldehyde OPA method.
- Add 2 uL of assay buffer (no dye, no sample) to the lower pedestal. Lower the arm and click F3 or the Blank button. When the measurement is complete, lift the arm and use a dry laboratory wipe to blot the buffer from both the bottom and upper measurement surfaces. Use a fresh aliquot of buffer to verify a proper baseline.
- Under Measurement type, click on the Standards tab. Highlight the Reference standard.
- Mix the reference solution (assay buffer and dye, no sample) briefly and transfer 2 uL of the solution onto the lower pedestal. Lower the arm and click F1 or the Measure button. A pop up window will ask for confirmation of the units. (Recommended ng/mL or pg/uL)
- Measure up to 5 replicates of the reference solution using a fresh 2 uL aliquot for each measurement.
- Select Standard 1 to enter a value. Enter values for up to 7 standards.
- Mix the standard solution briefly and transfer 2 uL onto the lower pedestal. Lower the arm and click F1 or the Measure button. Measure up to 5 replicates of each standard using a fresh 2 uL aliquot for each measurement.
- Once the standard curve is completed, select the Standard Curve Type (Interpolation, Linear, 2° polynomial, 3° polynomial) that best fits the standards data set.
- Click on the Sample tab under Measurement Type, and enter the unknown samples' respective ID information. If a dilution of the unknown sample was made, enter the dilution factor in the box below the sample ID window.
- Add 2 ul of the sample and use the F1 key or click the Measure button to initiate the measurement cycle. Use a fresh aliquot of sample for each measurement.

### Low Range (1 ug/ml–50 ug/ml) Dilution Series

BSA (ug/ml)	Ave RFU (n=5)	St dev	%CV
0	199	8.4	21
0.8	380	12	14
1.6	630	35	29
3.1	1305	19	6
6.3	2462	50	6.7
12.5	4858	31	1.9
25	8523	32	1
50	14863	79	1.4

### High Range (15 ug/ml–1000 ug/ml) Dilution Series

BSA (ug/ml)	Ave RFU (n=5)	St dev	%CV
0	199	56	28
16	16	75	20
31	31	48	7.7
63	63	22	1.7
125	125	162	6.7
250	250	150	3
500	500	79	0.9
1000	1000	256	1.7