

1 Abstract

Cytokine levels are measured in many different experimental models, with samples obtained in a variety of biological matrices such as cell culture supernatant, serum, plasma, saliva, cerebrospinal fluid, and tissue homogenate. The most common method of cytokine analysis is enzyme-linked immunosorbent assay (ELISA). However, ELISA often requires high sample volume (at least 50 µL), which is difficult to obtain when working with small animals such as mice. Also, the ELISA method has a limited dynamic range and low throughput due to its numerous wash steps. Considering the limitations of ELISA technology, a faster approach was developed for cytokine analysis in complex biological samples. This novel analytical approach is a chemiluminescent homogeneous bead-based immunoassay that is performed without wash steps. The cytokine of interest is sandwiched between two highly specific antibodies bound to the beads. In presence of the analyte, the beads come into proximity resulting in an emission of light upon laser excitation. The amount of light emitted is directly proportional to the amount of analyte present. Sample volumes as low as 5 µL can be utilized in this sensitive detection platform. In addition, upper quantitation limits can extend to much higher concentrations than achievable using traditional ELISA assay. For example, mouse IL6 can be detected from 1 pg/mL to 100,000 pg/mL without dilution, whereas a comparable ELISA has a range of 8 pg/mL to 500 pg/mL. The application of this immunoassay to different types of biological samples was verified by performing spike recovery experiments, and also by comparison of the levels measured to the levels reported in the literature. Overall, this assay technology will aid researchers in evaluating cytokines from many different biological sources that, in some cases, have been problematic to measure in the past.

2 Materials and Methods

Mouse and rat IL6, TNFα and MCP-1 detection:

AlphaLISA® kits available from PerkinElmer: IL6 (mouse) (AL504), TNFα (mouse) (AL505) and CCL2/MCP-1 (mouse/rat) (AL509). Assay buffer is AlphaLISA Immunoassay Buffer (AL000C) for mouse IL6 (mIL6) and mouse TNFα (mTNFα), and NaCl Buffer (AL007C) for mouse MCP-1 (mMCP-1). All calibration curves were done using mouse analyte supplied in the kit. All 3 kits similarly detect mouse and rat cytokines.

Mouse/rat sample fluids:

All mouse/rat sample fluids were supplied as pools of individuals, non-medicated, non-immunized, from Bioreclamation LLC:

- Mouse bronchial lavage fluid (BALF): Strain CD-1, cat# MSE-BRONLEV
- Mouse serum: Strain BALB/C, cat# MSESRLM-BALB
- Rat amniotic fluid (AF): Strain Sprague Dawley, cat# RATAMNFL
- Rat cerebrospinal fluid (CSF): Strain Sprague Dawley, cat# RATCSF

Diluents tested:

AlphaLISA HiBlock Buffer: PerkinElmer, cat# AL004C.

Foetal bovine serum Premium (FBS): Wisent, cat# 080150.

Fluids tested as diluents were supplied as pools of individuals, non-medicated, non-immunized, from Bioreclamation LLC:

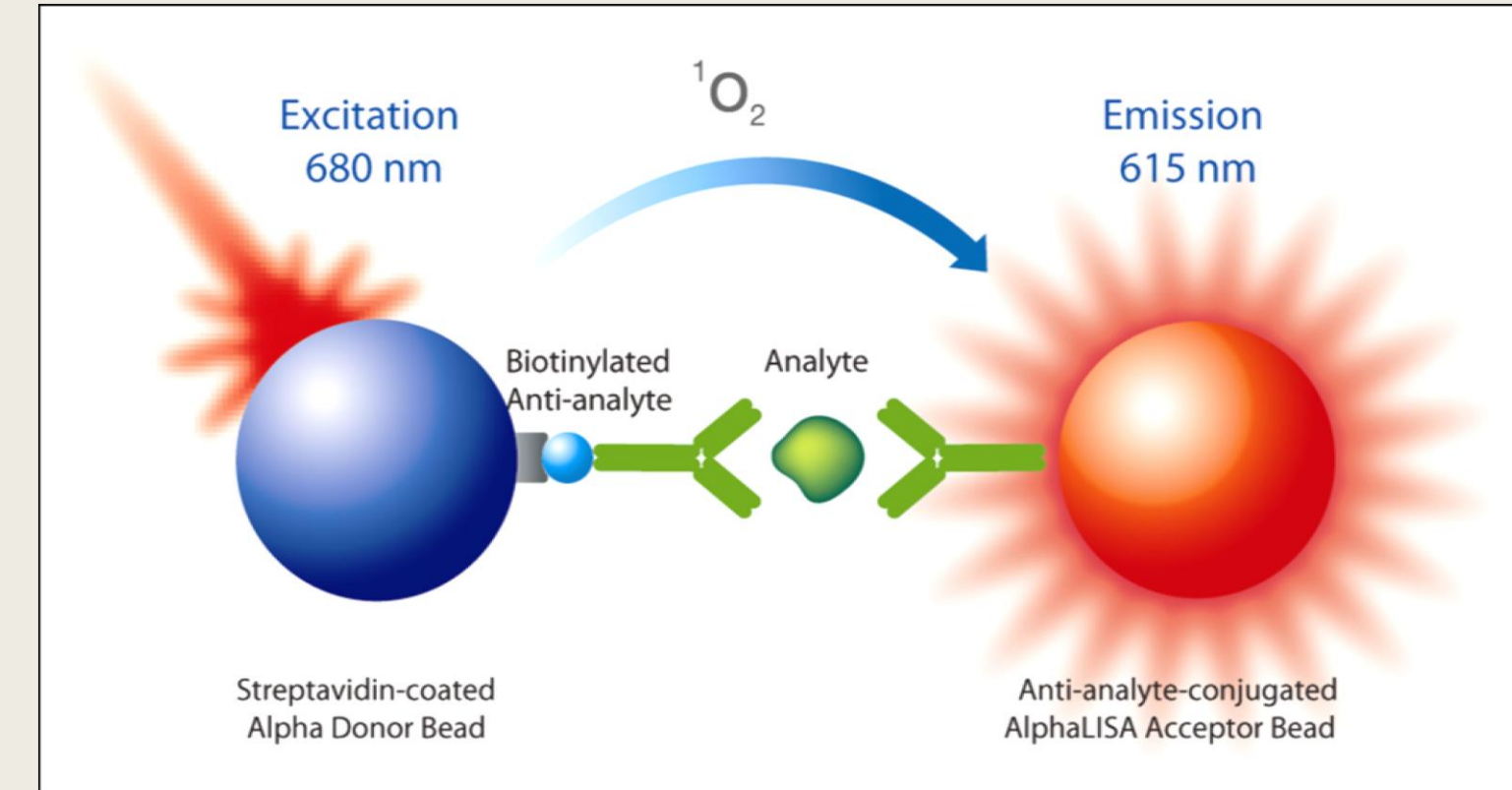
- Beagle bronchial lavage fluid (BALF): cat# BGL-BROLAV
- Beagle cerebrospinal fluid (CSF): cat# BGLCSF
- Rabbit amniotic fluid (AF): Strain New Zealand White, cat# RABAMNFL

Fluids pre-treatment and storage:

Prior to being used in AlphaLISA assays: FBS was triple filtered at 0.1µm by Wisent; mouse serum was filtered at 0.2µm by Bioreclamation LLC; BALF and AF fluids were centrifuged at 650xg for 15 min at 4°C to remove cells and cell debris, and supernatants were used in assays; CSF fluids were not pre-treated.

Storage: Mouse serum and FBS were kept at -20°C. All other fluids (BALF, CSF, AF) were kept at -80°C.

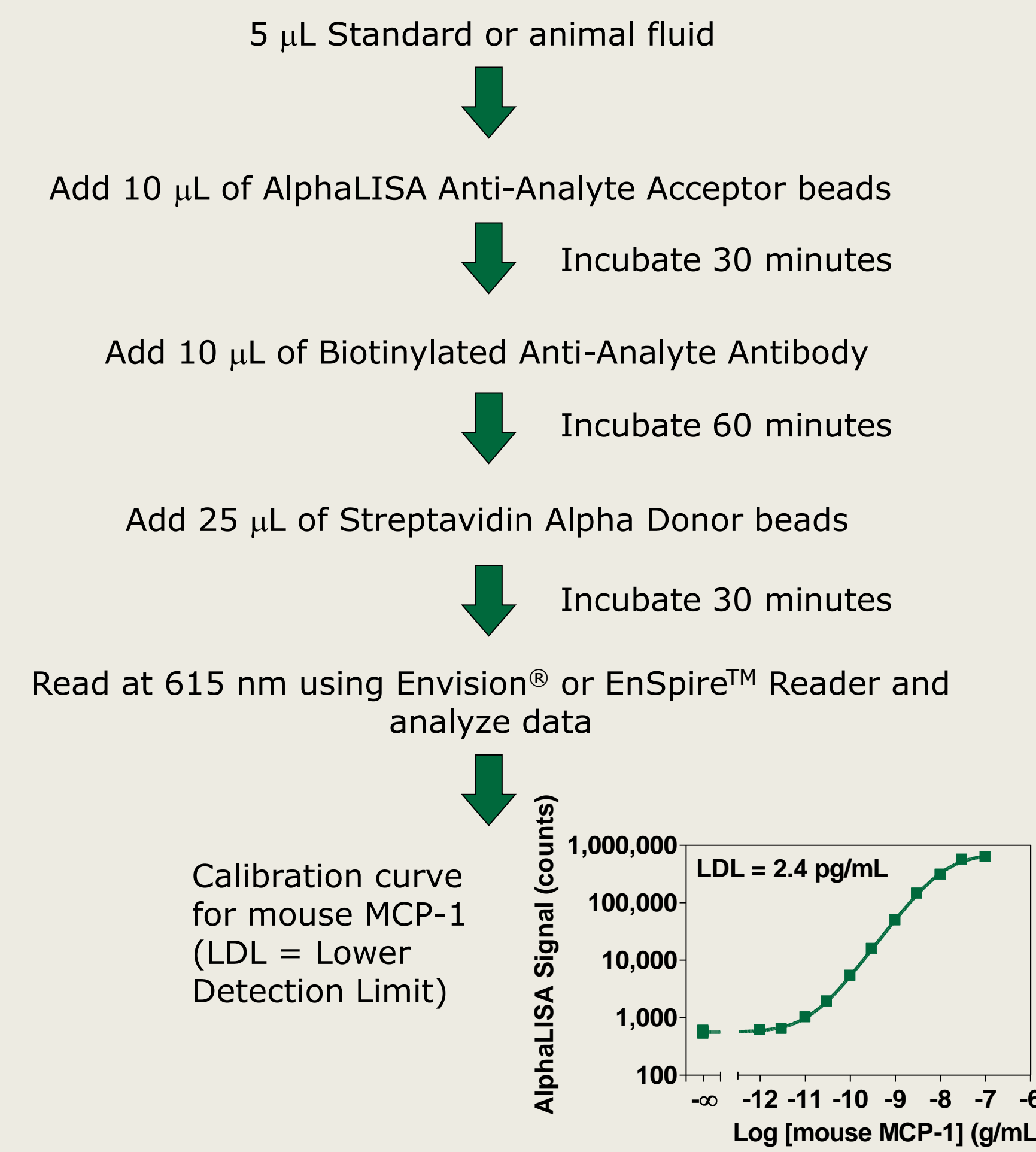
3 Alpha Technology Assay Principle



The biotinylated anti-analyte antibody binds to the Streptavidin-coated Donor beads while another anti-analyte antibody is conjugated to AlphaLISA Acceptor beads. In the presence of the analyte, the beads come into close proximity. The excitation of the Donor beads provokes the release of singlet oxygen molecules that triggers a cascade of energy transfer in the Acceptor beads resulting in a sharp peak of light emission at 615 nm.

4 Immunoassays – quick and simple

Assay protocol with 3 incubation steps (No wash steps)



5 Fluids and cytokines chosen

Cytokines and mouse/rat fluids were selected according to availability of mouse/rat PerkinElmer AlphaLISA kits, as well as availability and cost of fluids from suppliers.

Expected physiological range:

| Cytokine | Physiological range found in literature (pg/mL) | | | |
|----------|---|------------|-----------|---------|
| | Mouse serum | Mouse BALF | Rat CSF | Rat AF |
| IL6 | 5 - 70 | 20 | Not found | 10 - 60 |
| TNFα | 5 - 100 | 30-90 | Not found | 10 - 60 |
| MCP-1 | 30 - 1500 | 40 | Not found | 1500 |

6 Experimental approach

Step 1: Perform linearity experiment with the following diluents for all 3 cytokines:

| Tested diluent | Spiked sample fluid (3ng/mL or 10ng/mL mouse cytokine) | | | |
|---|--|------------|---------|--------|
| | Mouse serum | Mouse BALF | Rat CSF | Rat AF |
| Assay buffer from kit | X | X | X | X |
| HiBlock buffer | X | X | X | X |
| PBS-0.1%BSA | X | X | X | X |
| PBS-0.01%BSA | X | X | X | X |
| Same fluid as sample fluid, but other species | X | X | X | X |
| FBS* | X | | | |

*For serum samples, FBS was the only diluent tested, as FBS has previously been shown to be the best diluent for all serum samples (data not shown).

Step 2: Diluent selection:

With step 1 results, select the two best diluents based on:

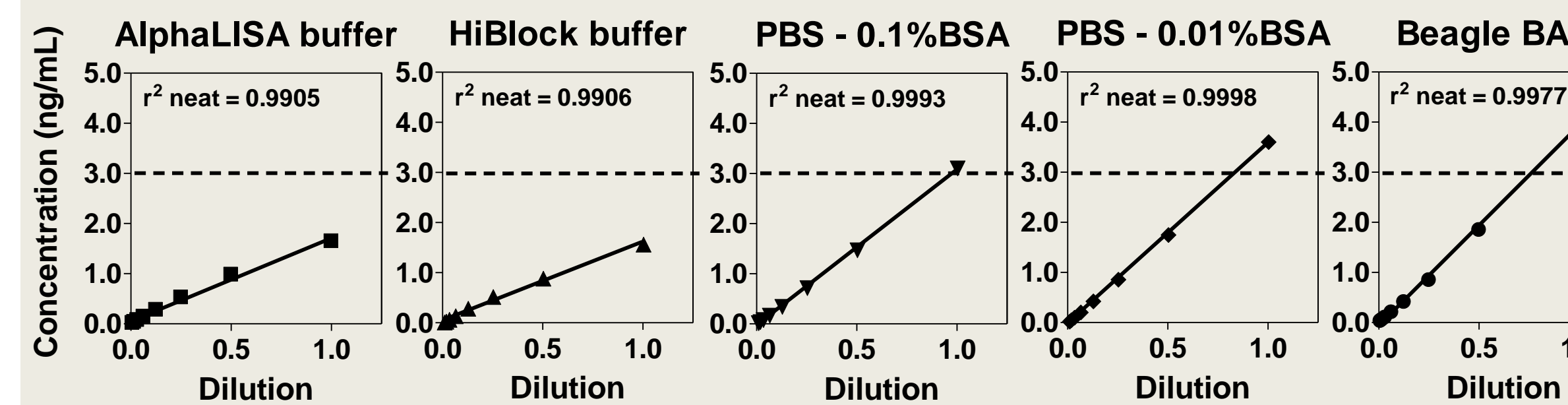
- Observed concentration X dilution factor (DF) as close as possible to spiked value, for all dilution points
- Recoveries between 70-130% for all dilution points

Step 3: Evaluate spike and recovery value:

- Spike neat and 1/2 diluted mouse/rat fluid with mouse cytokine
- Use spiked diluent as spike control
- Select 3-4 spikes and include a "no spike" control (to measure physiological level)
- Recoveries between 70-130% are targeted
- Evaluate other diluents if good recovery is not obtained

7 Example (no dilution required)

Step 1: Linearity of dilutions in 5 diluents for spiked mouse BALF (3ng/mL mTNFα)



| Dilution factor (DF) | AlphaLISA buffer | HiBlock buffer | PBS-0.1%BSA | PBS-0.01%BSA | Beagle BALF |
|----------------------|------------------------|------------------------|------------------------|------------------------|------------------------|
| | Recovery from neat (%) | Recovery from neat (%) | Recovery from neat (%) | Recovery from neat (%) | Recovery from neat (%) |
| 1 | 100 | 100 | 100 | 100 | 100 |
| 2 | 119 | 113 | 95 | 97 | 92 |
| 4 | 128 | 135 | 93 | 96 | 85 |
| 8 | 137 | 150 | 90 | 96 | 82 |
| 16 | 139 | 146 | 92 | 95 | 83 |
| 32 | 149 | 145 | 94 | 99 | 83 |

Step 2:

- Diluents not chosen: Linearity not good from neat
- Choose 2 good diluents: PBS-0.1%BSA and Beagle BALF
- Good linearity from neat
- Observed concentration X DF close to spiked value (3ng/mL) for all dilutions

Step 3: Chosen diluent showing the best spike and recovery results

| Spike (pg/mL) | Diluent: PBS-0.1%BSA | | |
|---------------|------------------------|------------------------|----------------|
| | Spiked diluent control | Spiked neat Mouse BALF | Recovery (%)** |
| No spike | 0.0 | 17.1 | NA |
| 10 | 12.0 | 9.2 | 76 |
| 30 | 35.2 | 32.3 | 92 |
| 300 | 300.2 | 292.6 | 97 |
| 3000 | 3141.2 | 2952.1 | 94 |

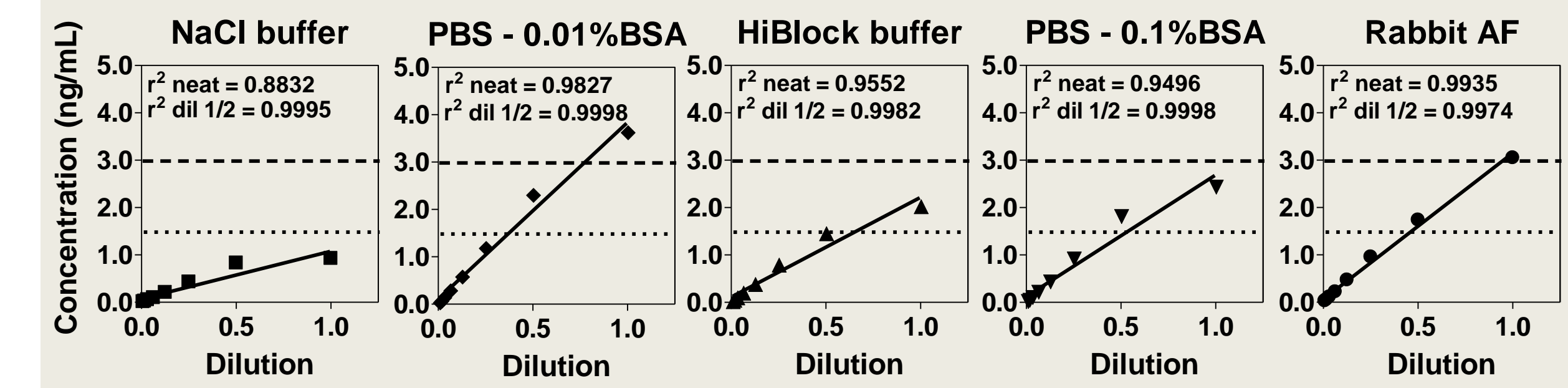
Excellent linearity of dilutions from neat (Step 1) and excellent recovery for all 4 spikes tested. Measured TNFα physiological level in mouse BALF (No spike) = 17.1 pg/mL

* Concentration for 10-3000 pg/mL spikes = measured concentration – no spike value.

** Recovery (%) = recovery compared to spiked diluent control

8 Example (1/2 dilution required)

Step 1: Linearity of dilutions in 5 diluents for spiked rat AF (3ng/mL mMCP-1)



| Dilution factor (DF) | NaCl buffer | PBS-0.01%BSA | HiBlock buffer | PBS-0.1%BSA | Rabbit AF |
|----------------------|------------------------|------------------------|------------------------|------------------------|------------------------|
| | Recovery from neat (%) | Recovery from neat (%) | Recovery from neat (%) | Recovery from neat (%) | Recovery from neat (%) |
| 1 | 100 | 100 | 100 | 100 | 100 |
| 2 | 179 | 100 | 144 | 149 | 114 |
| 4 | 187 | 104 | 157 | 152 | 126 |
| 8 | 180 | 100 | 129 | 101 | 145 |
| 16 | 170 | 95 | 129 | 101 | 163 |
| 32 | 192 | 107 | 129 | 101 | 169 |

Step 2:

- Diluents not chosen: observed concentration X DF not close to spiked value (3ng/mL) (too low or too high).
- Choose 2 good diluents: PBS-0.1%BSA and rabbit AF
- Excellent linearity from 1/2 dilution
- Observed concentration X DF close to spiked value (3ng/mL) for all dilutions.

Step 3: Chosen diluent showing the best spike and recovery results:

| Spike (pg/mL) | Diluent: PBS-0.1%BSA | | | |
|---------------|----------------------|--------------------|-------------------|---|
| | Spiked diluent | Spiked neat Rat AF | Spiked 1/2 Rat AF | 1/2 dilution in diluent required: excellent recovery for all 3 spikes tested in 1/2 rat AF, but not in neat rat AF. Measured MCP-1 physiological level in rat AF (No spike) = 575.6 pg/mL X 2 = 1151pg/mL |
| No spike | 0.0 | 910.5 | NA | 575.6 |
| 300 | 267.3 | 195.2 | 73 | 317.2 |
| 1000 | 890.5 | 442.6 | 50 | 750.4 |
| 3000 | 2627.6 | 1542.0 | 59 | 2227.9 |

* Concentration for 300-3000 pg/mL spikes = measured concentration – no spike value.

** Recovery (%) = recovery compared to spiked diluent control

9 Summary table

| Cytokine | Sample fluid | Best diluent | LDL in diluent (pg/mL) | Dilution of sample fluid required in diluent | Physiological level measured (pg/mL) | Spike range in sample fluid giving good recovery (pg/mL) |
|----------|--------------|--------------|------------------------|--|--------------------------------------|--|
| IL6 | Mouse serum | FBS | 2.1 | 1/2 | 33 | 30 - 10000 |
| | Mouse BALF | PBS-0.1%BSA | 2.4 | Neat | 22 | 10 - 3000 |
| | Rat CSF | Beagle CSF | 3.4 | Neat | ND | 10 - 3000 |
| TNFα | Mouse serum | FBS | 1.6 | 1/2 | 55 | 30 - 3000 |
| | Mouse BALF | PBS-0.1%BSA | 1.6 | Neat | 17 | 10 - 3000 |
| | Rat CSF | PBS-0.1%BSA | 1.4 | Neat | ND | 3 - 3000 |
| MCP-1 | Mouse serum | FBS | 5.4 | 1/2 | 444 | 300 - 3000 |
| | Mouse BALF | Beagle BALF | 5.7 | Neat | 22 | 10 - 3000 |
| | Rat CSF | PBS-0.1%BSA | 2.9 | Neat | 501 | 100 - 3000 |
| | Rat AF | PBS-0.1%BSA | 2.4 | 1/2 | 1151 | 300 - 3000 |

10 Conclusion

- Sensitive bead-based immunoassays were developed to detect IL6, TNFα and MCP-1 from rodent serum, BALF, CSF and AF. Good recoveries were obtained from fluids spiked as low as 3 pg/mL.
- This study highlights that the appropriate diluent is not easy to predict. In some cases where a perfect match of diluent and fluid is not found, sample dilution is required.
- These no wash rapid assays require only 5µL sample volumes. This represents a substantial improvement over ELISA-type assays in terms of speed of execution and sample consumption.
- Moreover, results presented demonstrate that measured physiological levels are also in agreement with the literature.