

Kayla J. Smith<sup>1</sup>, Bruce A. Sherf<sup>1</sup>, Francesca Ferguson<sup>2</sup>, Heather E. Preisendanz<sup>3</sup> and John P. Vanden Heuvel<sup>1,4</sup>

<sup>1</sup>INDIGO Biosciences, Inc., State College, PA 16801, USA, <sup>2</sup>Department of Entomology, The Pennsylvania State University, University Park, PA 16802, USA, <sup>3</sup>Department of Agricultural and Biological Engineering, The Pennsylvania State University, University Park, PA 16802, USA, <sup>4</sup>Department of Veterinary and Biomedical Sciences, The Pennsylvania State University, University Park, PA 16802, USA

#18948

## Abstract

Current protocols for water safety evaluation primarily focus on targeted chemical analysis, but this approach does not alert us to the presence of hazardous bioactivities from non-target chemical contaminants. Importantly, the cumulative bioactivities from low concentrations of numerous mixed chemical pollutants cannot be predicted using targeted analytical approaches. As a complement to analytical chemistry approaches, mammalian cell-based assays have been developed to provide a sensitive assessment of potential adverse effects from chemical contaminants in water such as endocrine-disruption (ED). Herein we demonstrate the utility of new cell-based luciferase reporter assays encompassing chemical-sensing receptors that are directly relevant to environmental biomonitoring applications. These include the polycyclic aromatic hydrocarbon (PAH) sensing Aryl Hydrocarbon Receptor (AhR), and the ED targets Androgen Receptor (AR), Estrogen Receptor (ER), Glucocorticoid Receptor (GR), and the Mineralocorticoid Receptor (MR). Environmentally relevant reference ligands for each receptor were tested in the assays, evaluated for sensitivity, and compared to current or suggested monitoring trigger levels (MTLs). Surface water samples collected from nine sites in the Spruce Creek watershed in Central Pennsylvania were tested in these assays. The assay data were compared to targeted chemical analyses of the sites to determine if the assays were able to detect biological activities that the chemical data would not have predicted. The results demonstrate that the limit of detection (LOD) for each assay is consistent with proposed MTLs presented in the literature. All samples exhibited detectable agonist activities for AhR, AR, ER, and MR, but minimal to no activity for GR. The comparison between bioassay and chemical analysis data demonstrated that using both methods together provide a better representation of the quality of the water samples, and that chemical analysis alone cannot predict the accumulated bioactivities revealed by the bioassays. These results demonstrate the utility and sensitivity of function-based assays for generating comprehensive water quality assessments. Bioassays provide reliable, actionable data that inform the decision-making process as to the need and extent of remediation required for a given water source.

## Materials and Methods

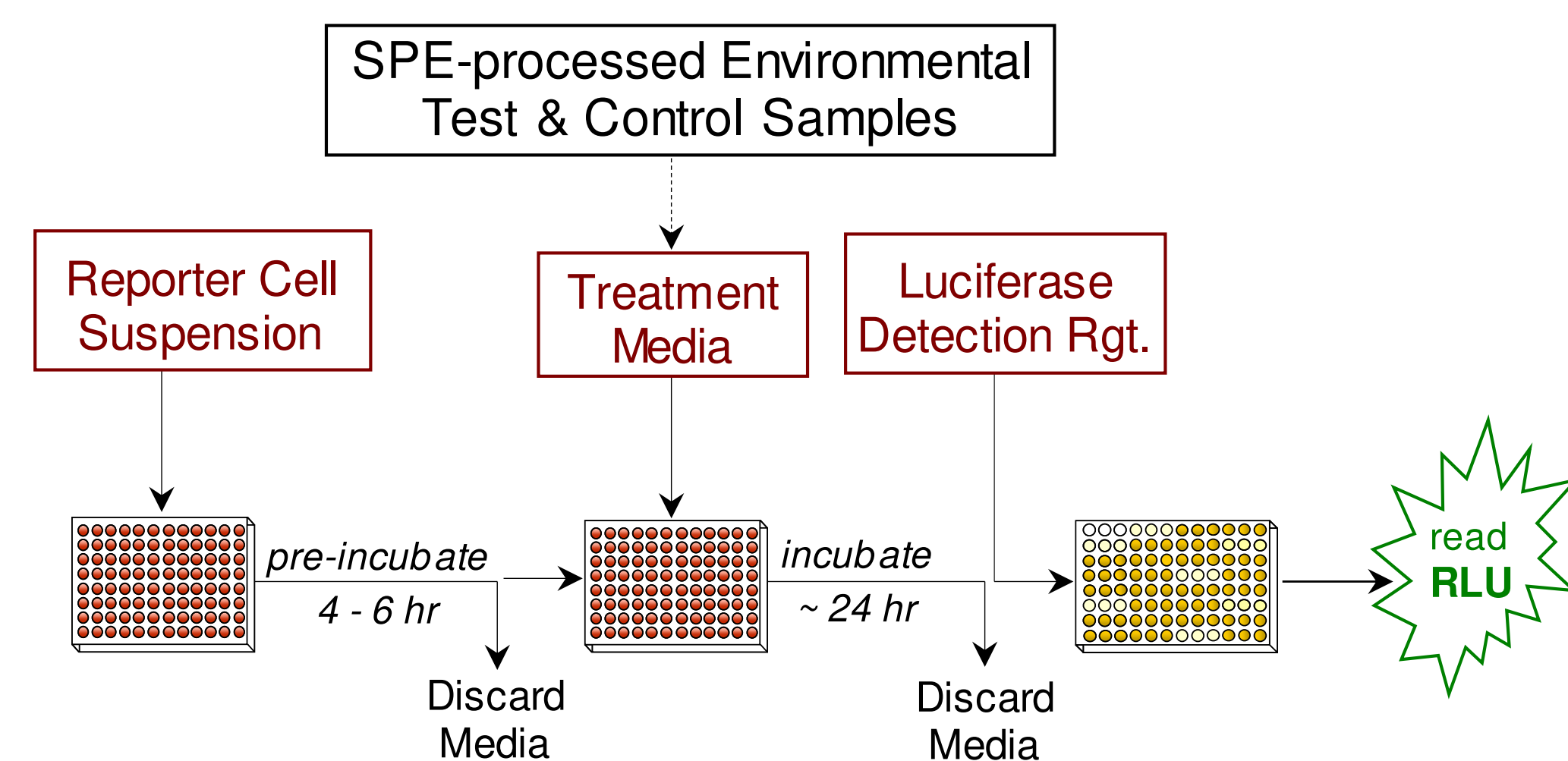
**Sample collection and processing:** Water samples were collected on one day in August 2021 from nine sites in the Spruce Creek watershed in central Pennsylvania, plus one field blank. These groundwater collection sites represent mixed use water sources (forested 59%, agriculture 35%, developed 6%). Water samples were processed before bioassay analyses *via* solid-phase extraction (SPE) using 200 mg Oasis® HLB sorbents (Waters Corp.) and elution solvent mixes, as described by Vanden Heuvel, *et. al.* (Science of the Total Environment, manuscript in preparation).

**Bioassay setup:** Luciferase reporter assays were developed and optimized in mammalian cell lines for the following receptors: AhR, AR, ER, GR, and MR. Figure 1 depicts the setup and workflow of the five environmental bioassays.

**Assay data analyses:** For each assay a standard curve was generated using an environmentally relevant reference analyte. Each assay's limit of detection (LOD) was determined using an unpaired, two-tailed t-test to assess changes in RLU values between the vehicle-treated cells and the low-concentration reference agonist-treated cells. Significance was set at  $p \leq 0.01$ . Field Blank RLU values were used for background-subtraction from the Test Sample RLU values. For each Test Sample, values of Percent Relative Activation were calculated by normalizing RLU's to the maximal RLU value achieved by each receptor's reference agonist (= 100%). Respective standard curves were then used to interpolate bio-equivalent (BEQ) concentrations of each test sample.

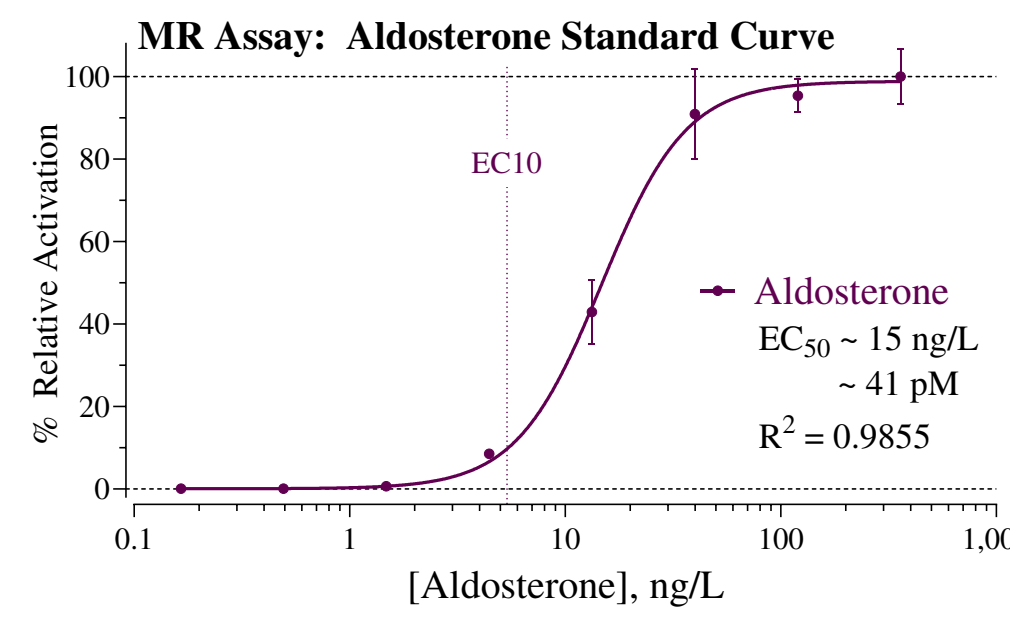
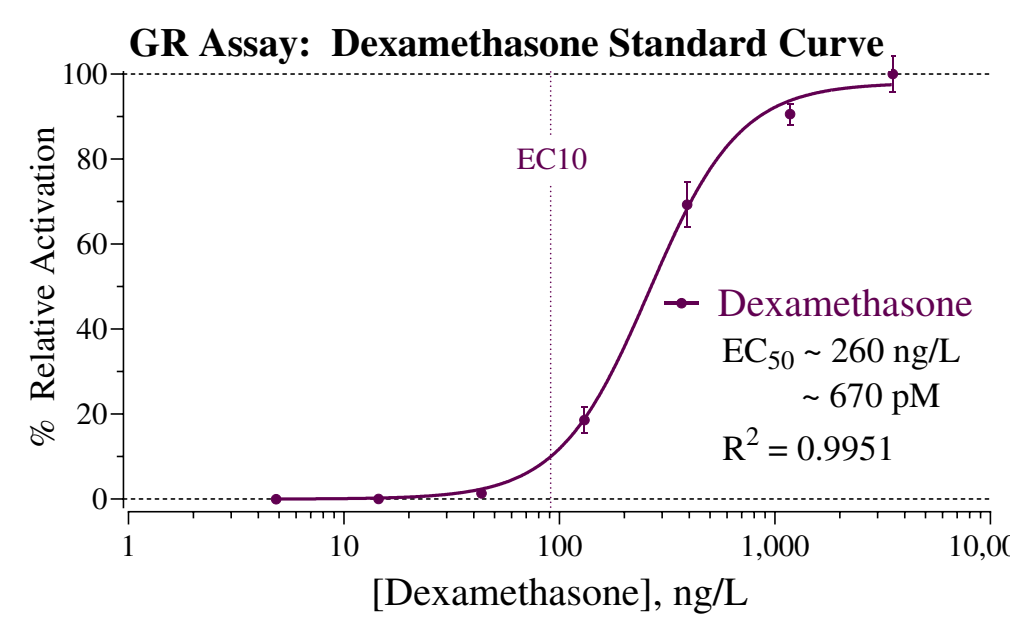
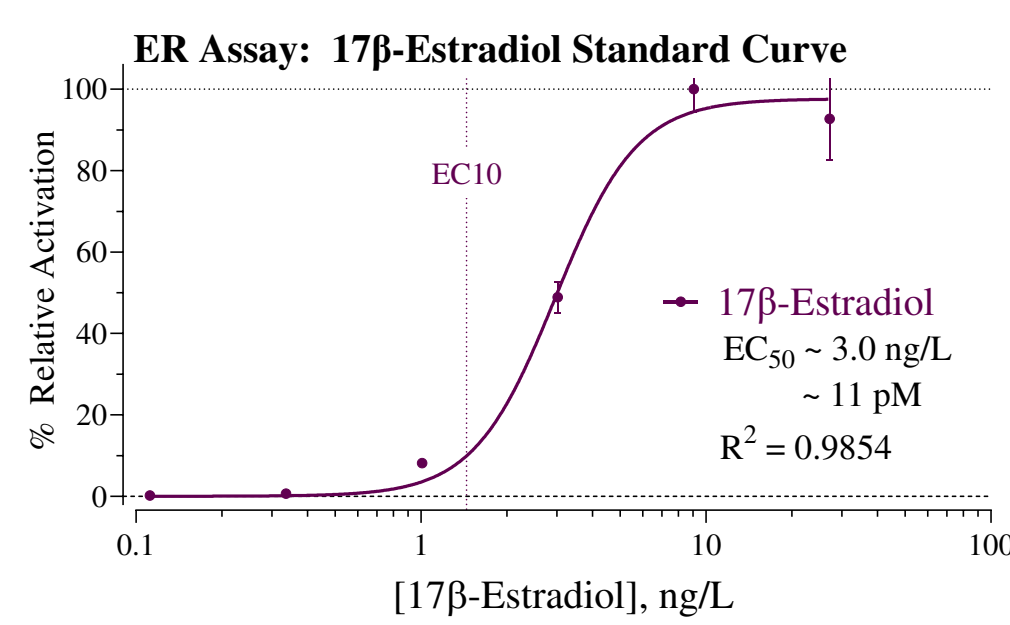
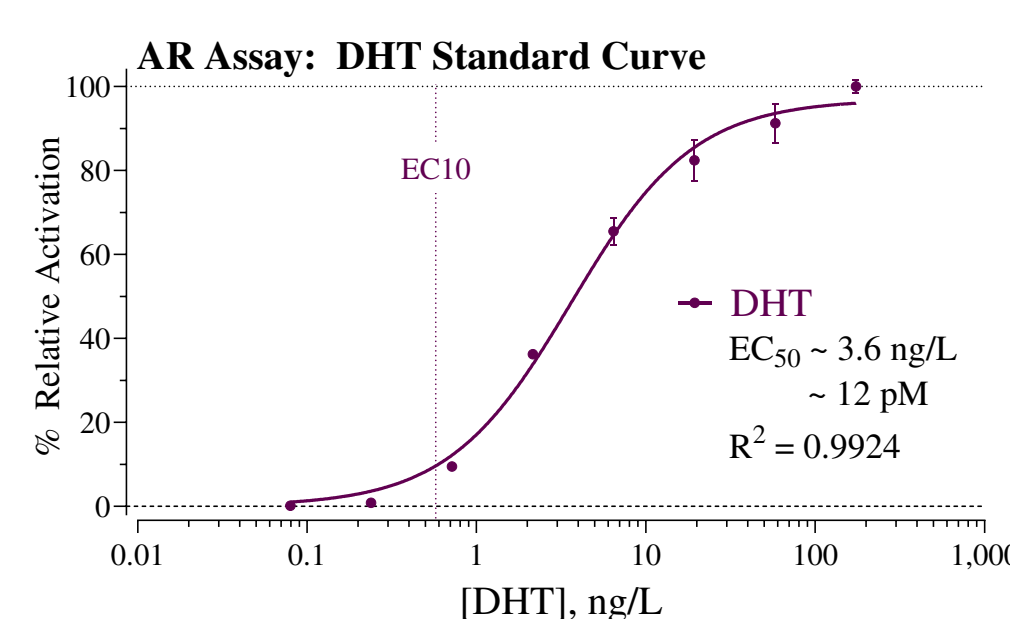
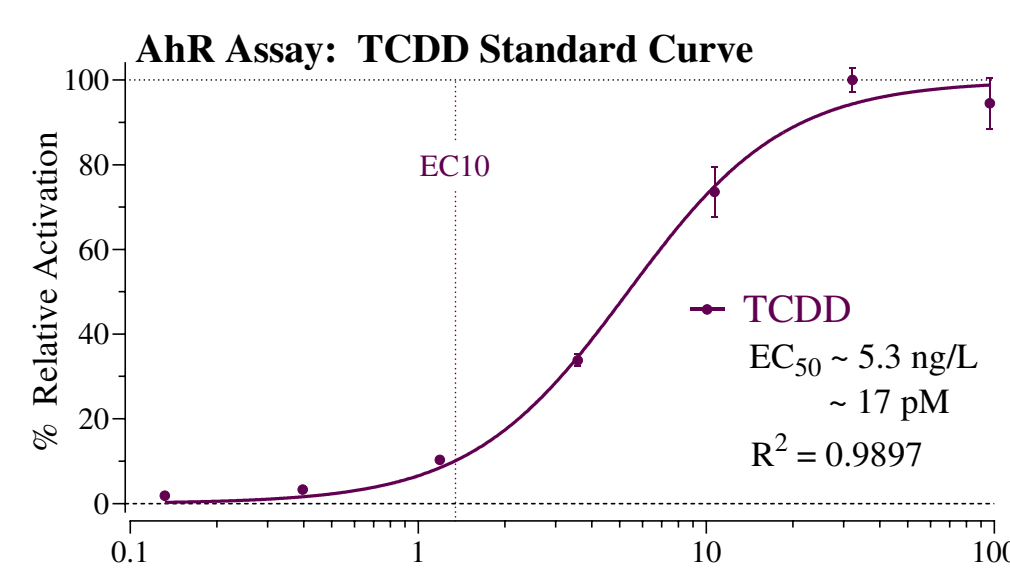
**Chemical analyses:** SPE-processed water samples were also analyzed for chemical contaminants including pharmaceuticals, pesticides, and personal care product components. Analyses were conducted using high-resolution accurate mass (HRAM) Q Exactive mass spectrometer interfaced with an ICS-5000+ chromatography system *via* a heated electrospray injection source, as described in Vanden Heuvel, *et. al.* (Science of the Total Environment, manuscript in preparation).

**References:** \*California State Water Resources Control Board, Water Quality Control Policy for Recycled Water, Adopted December 11, 2018. #Values summarized in Neale, *et. al.*, (2023) Effect-Based Trigger Values Are Essential for the Uptake of Effect-Based Methods in Water Safety Planning. Environmental Toxicology and Chemistry. 42: 714-726. Values depicted are those for Human EBT-BEQ (drinking and recycled water for indirect potable reuse).

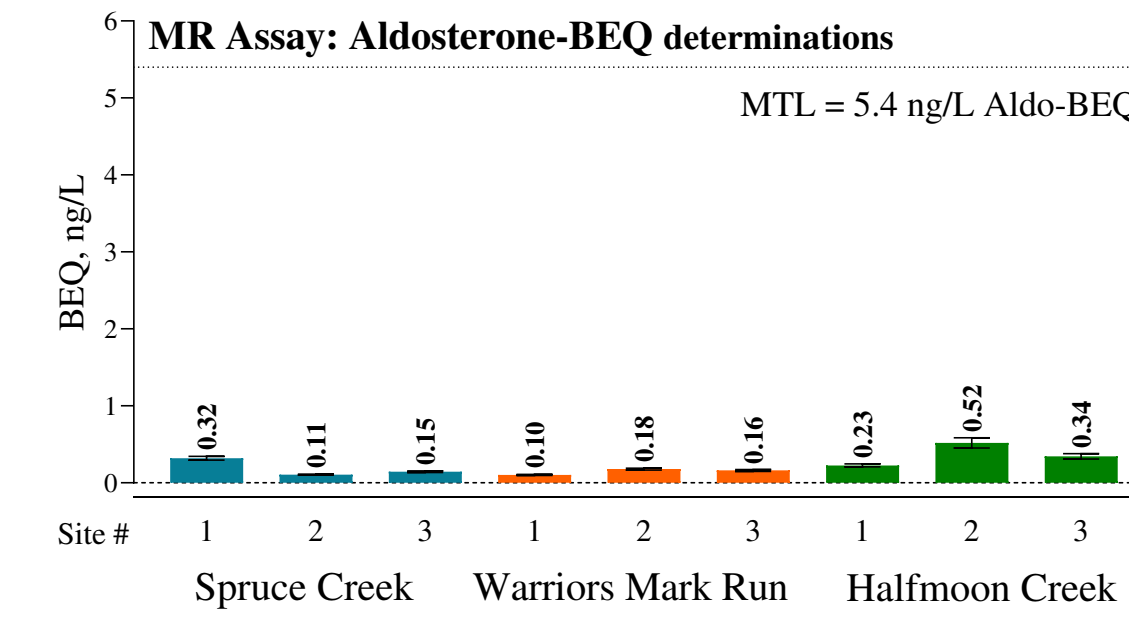
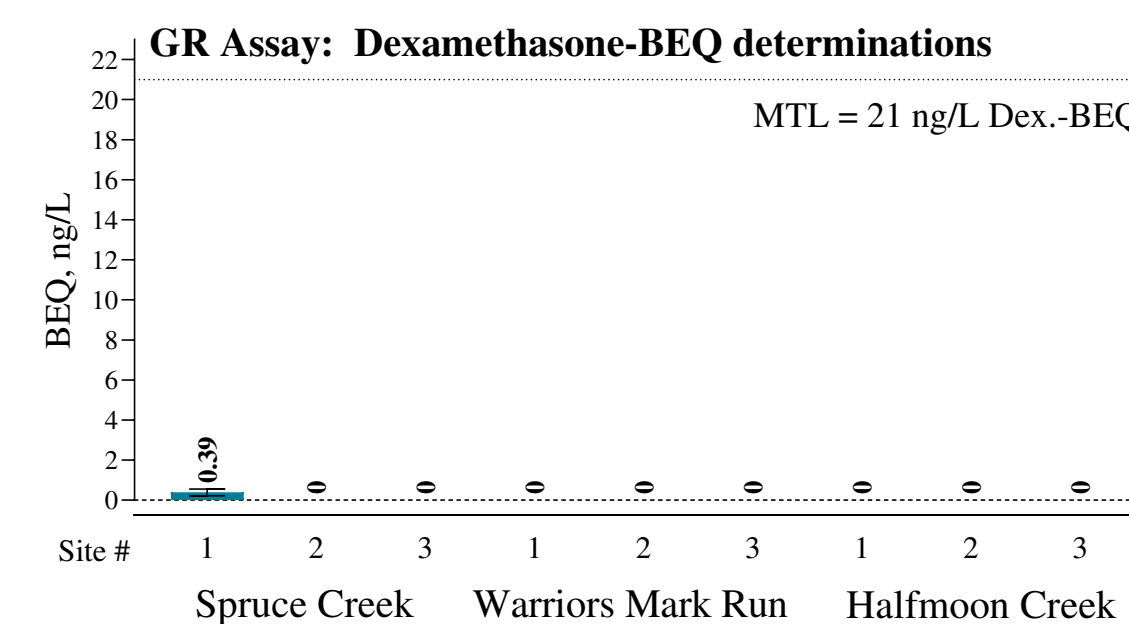
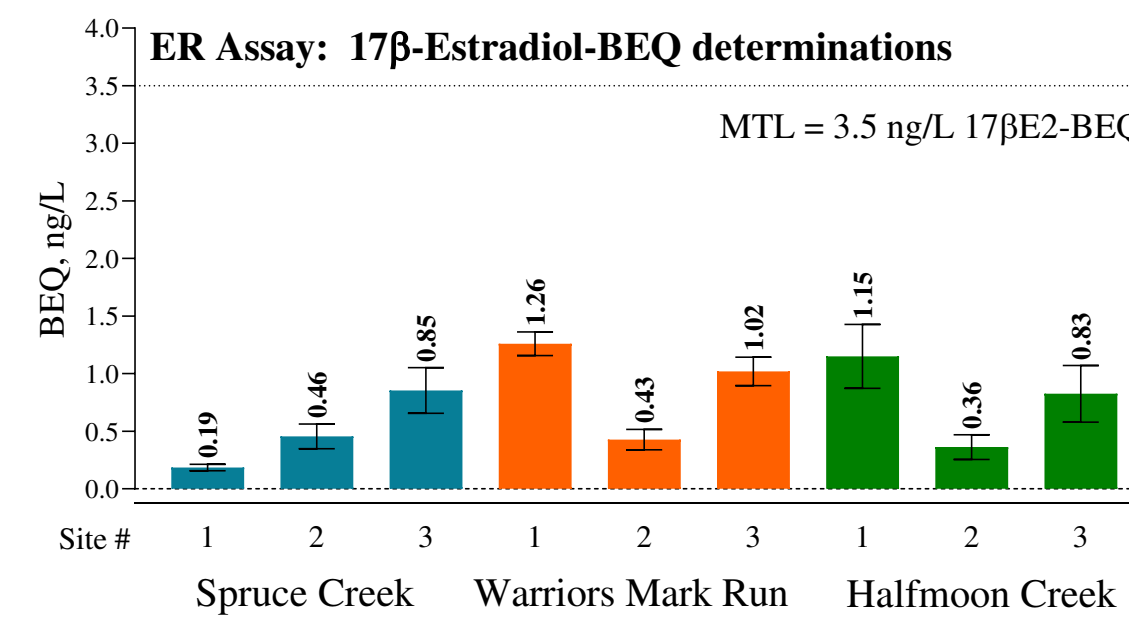
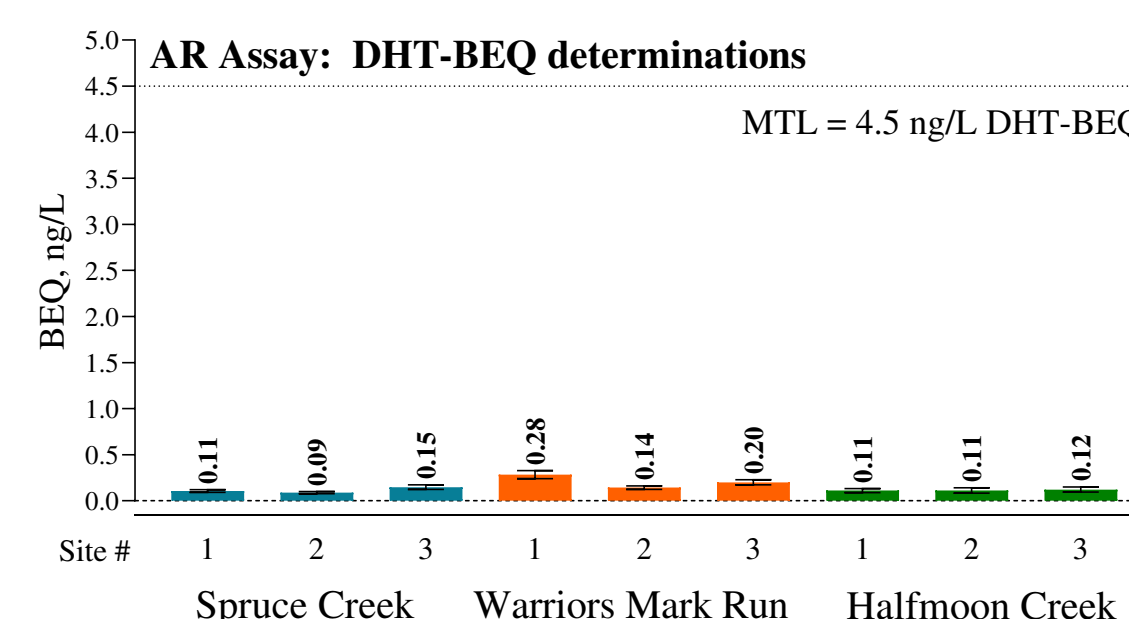
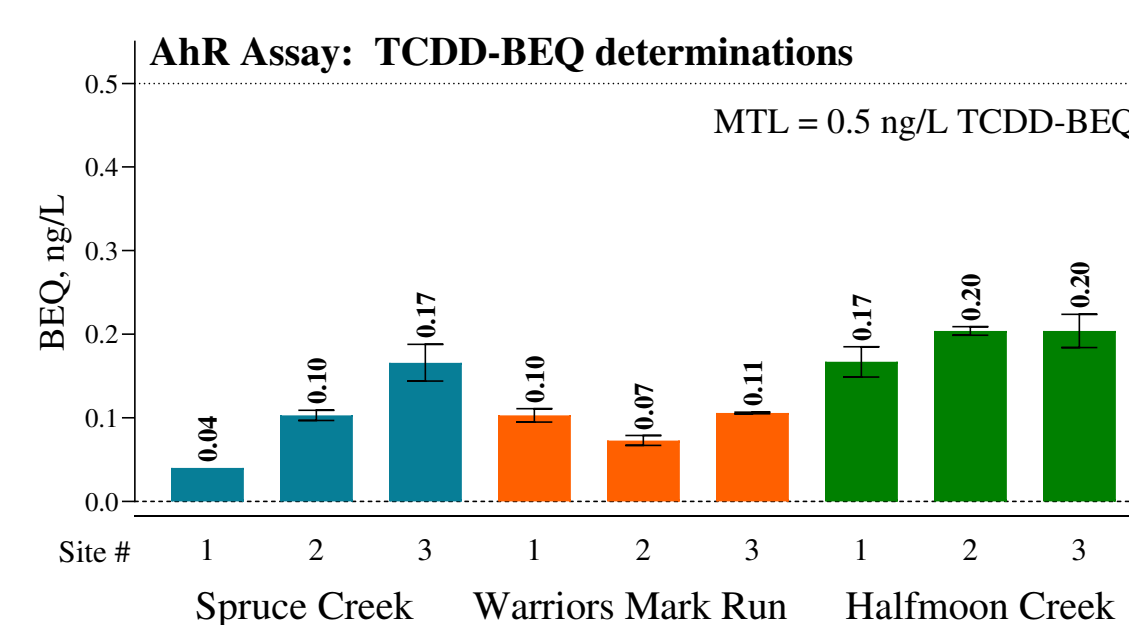


**Figure 1: Bioassay setup and workflow.** Assay protocols were performed as prescribed by INDIGO Biosciences (State College, PA, USA)

**Figure 2**



**Figure 3**



## Results and Conclusions

**Figure 2: Standard Curves for the AhR, AR, ER, GR and MR bioassays.** Dose-response curves are generated using environmentally-relevant reference agonists.

**Figure 3: BEQ determinations for water samples using the AhR, AR, ER, GR, and MR bioassays.** The Standard Curves depicted in Figure 2 were used to extrapolate respective BEQ values for each field sample tested. MTL values for each reference are depicted and summarized in Table 1. For the MR assay the EC<sub>10</sub> value for Aldosterone is used as the MTL. Field sites that yield test sample BEQ values exceeding the MTL value require further investigation and potential remediation.

Assay	Reference Compound	EC <sub>50</sub>	EC <sub>10</sub>	LOD	MTL
AhR	TCDD	5.3	1.3	0.40	0.5*
AR	DHT	3.6	0.58	0.24	4.5 - 32#
ER	17β-Estradiol	3.0	1.4	1.0	3.5*
GR	Dexamethasone	260	90	15	21 - 150#
MR	Aldosterone	15	5.4	0.49	na

**Table 1: Summary of Environmental Bioassay Metrics.**

Assay metrics for the AhR, AR, ER, GR and MR assays are depicted using the reference agonists indicated.

na = not available  
TCDD = 2,3,7,8-Tetrachlorodibenzo-*p*-dioxin  
DHT = dihydrotestosterone

Chemical name	Field Blank	Spruce Creek			Warriors Mark Run			Halfmoon Creek			Class	Type	
		1	2	3	1	2	3	1	2	3			
Acetaminophen	0.0717	0.0059	0.0027	0.0010	0.0023	0.0066	0.0015	0.0031	0.0029	0.0008	Personal Care	NSAID	
Benzophenone-1	0.0025	0.0001	0.0001	0	0.0001	0	0	0	0.0002	0		Sunscreen	
Benzophenone-3	0.0106	0.0032	0.0001	0.0007	0	0.0001	0.0004	0.0001	0	0.0001		Sunscreen	
Caffeine	0.0007	0.0061	0.0213	0.0124	0.0056	0.0063	0.0070	0.0298	0.0031	0.0217	Personal Care	Stimulant	
DEET	0.0002	0.0004	0.0006	0.0013	0.0004	0.0011	0.0003	0.0017	0.0006	0.0016		Insect repellent	
Ibuprofen	0	0.192	0.956	0	0.142	0.077	0	1.24	0	0.191		NSAID	
Ketoprofen	0	0.0001	0	0.0009	0.0001	0	0.0001	0.0001	0.0001	0.0001	Personal Care	NSAID	
Naproxen	0	0.0009	0.0005	0.0007	0.0003	0.0006	0.0011	0.0003	0.0009	0.0013		NSAID	
Sucralose	0.0005	0	0.0004	0	0	0	0	0	0	0		Sweetener	
Theobromine	0.0005	0.0034	0.0114	0.0075	0.0038	0.0042	0.0076	0.022	0.0036	0.0159	Personal Care	Cosmetics	
Atrazine	0.0008	0.0691	0.0425	0.0357	0.0266	0.0248	0.0469	0.0203	0.0286	0.0213		Pesticide	Herbicide
Carbaryl	0	0.0036	0.0035	0.0041	0	0	0	0	0	0.0002			Carbamate insecticide
Chlorpyrifos	0.0146	0.0078	0.0040	0.0042	0.0024	0.0009	0.0012	0.0011	0	0.0007	Organophosphate insecticide		
Clothianidin	0	0.0043	0.0033	0.0046	0.0045	0.0046	0.0073	0.0025	0.0087	0.0042	Pesticide	Neonicotinoid insecticide	
Cyhalothrin	1.17	0	1.30	0	0	0	1.30	0	0	0		Pyrethroid insecticide	
Esfenvalerate	0.991	0	1.02	1.02	1.02	1.02	1.02	1.02	1.02	1.02		Pyrethroid insecticide	
Imidacloprid	0	0.0028	0.0052	0.0109	0.0019	0.0017	0	0.0091	0	0.0085	Pesticide	Neonicotinoid insecticide	
Malathion	0.0008	0.0006	0.0001	0	0	0	0	0	0	0		Organophosphate insecticide	
Permethrin	3.15	3.14	3.14	3.15	3.15	3.15	3.15	3.15	3.15	3.15		Pyrethroid insecticide	
Simazine	0.0002	0.0029	0.0026	0.0032	0.0020	0.0025	0.0015	0.0019	0.0024	0.0022	Pesticide	Herbicide	
Ampicillin	0	0	0	0.0025	0.0008	0	0	0.0075	0	0.0027		Pharmaceutical	Antibiotic
Atenolol-MRT	0	0.0115	0	0	0	0	0	0	0	0			Beta Block
Bupropion	0	0.0129	0.0021	0.0014	0.0010	0	0	0	0	0	Antianxiety		
Carbamazepine-A	0.0004	0.0001	0.0076	0.0035	0.0016	0.0013	0	0.0298	0	0.0195	Pharmaceutical	Anticonvulsant	
Chlortetracycline	0.0007	0.0021	0.0017	0.0016	0.0029	0.0014	0.0021	0.0018	0	0		Antibiotic	
Citalopram	0	0.0194	0.0001	0.0001	0.0003	0.0001	0.0006	0.0002	0.0011	0.0003		Antidepressant	
Clarithromycin	0.0003	0.0020	0	0	0	0	0	0	0	0	Pharmaceutical	Antibiotic	
Erythromycin	0	0.0002	0	0	0	0	0	0	0	0		Antibiotic	
Fluoxetine	0.632	0.631	0.633	0.634	0.634	0.634	0.634	0.634	0.634	0.634		Antidepressant	
Metformin	0.0001	0.0007	0.0001	0.0002	0.0002	0.0001	0.0001	0	0	0	Pharmaceutical	Antidiabetic	
Metoprolol	0.0001	0.0038	0.0018	0	0.0023	0.0030	0.0017	0.0037	0	0.0008		Blood pressure	
Ofloxacin	0	0.0239	0.0132	0.0017	0.0029	0.0014	0.0003	0.0009	0.0001	0		Antibiotic	
Oxytetracycline	0	0.0020	0.0013	0.0006	0.0016	0.0018	0.0033	0.0015	0.0015	0.0036	Pharmaceutical	Antibiotic	
Sulfadiazine	0	0.0003	0	0	0	0	0	0	0	0		Antibiotic	
Sulfadimethoxine	0	0.0001	0	0	0	0	0	0	0	0		Antibiotic	
Sulfamethazine	0	0	0	0	0	0	0.0004	0	0	0	Pharmaceutical	Antibiotic	
Sulfamethoxazole	0	0.0021	0.0005	0.0001	0	0	0.0001	0.0011	0	0.0007		Antibiotic	
Tetracycline	0	0.0012	0.0007	0	0	0	0	0	0	0		Antibiotic	
Trimethoprim	0.0001	0.0170	0.0027	0.0010	0.0009	0.0002	0.0004	0.0056	0.0003	0.0042	Pharmaceutical	Antibiotic	
Venlafaxine	0	0.0057	0.0024	0.0014	0.0009	0.0005	0.0009	0.0046	0.0003	0.0034		Antidepressant	

### Conclusions:

- Bioassays are sensitive tools for detecting synergistic activities of complex chemical mixtures in environmental water samples.
- Targeted chemical analysis alone cannot predict the activities observed in the bioassays.
- Combined data from cell-based bioassays and targeted chemical analysis provide a broader picture of water sample quality, leading to more effective decision making.

**Table 2: Chemical analyses of water test samples.**

Concentrations of chemicals detected in the water samples are presented as mg/L; values are reported as  $\leq 3$  significant figures.

Also analyzed but not detected in any sample:

- 4-t-Octylphenol
- β-cyfluthrin
- Bifenthrin
- Thiacloprid
- Thiamethoxam,
- Thiamethoxam-d3
- Amoxicillin