

# Phosphoproteomics reveals GITR agonism differentially regulates inflammatory vs. regulatory T cell subsets.

Andrew Davis<sup>1</sup>, Lori Cunningham<sup>1</sup>, Matthew Stokes<sup>1</sup>, Alissa Nelson<sup>1</sup>, Kathryn Abell<sup>1</sup>, Michael Lewis<sup>1</sup>, Roy Scialdone<sup>1</sup>, Than Bui<sup>1</sup>, Emily Alonzo<sup>1</sup>, Gordon Freeman<sup>2,3</sup>, Sarah R Klein<sup>1</sup>  
<sup>1</sup>Cell Signaling Technology, Inc., <sup>2</sup>Dana-Farber Cancer Institute, <sup>3</sup>Harvard Medical School

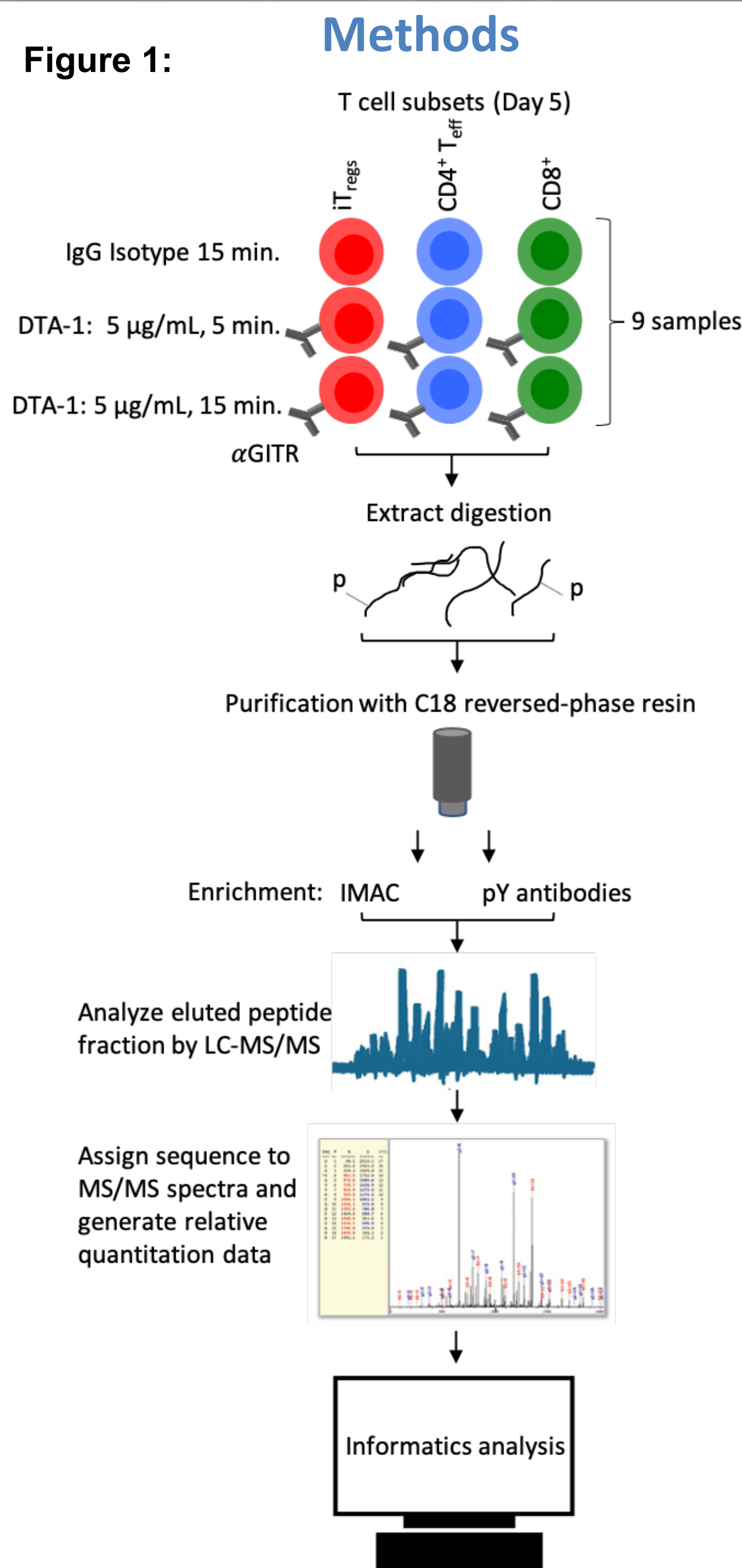
## INTRODUCTION

Triggering the co-stimulatory receptor glucocorticoid-induced TNFR-related protein (GITR) on T cells has been identified as a promising cancer immunotherapeutic strategy, yet the signaling events by which GITR agonism induces antitumor immunity are not well understood. GITR is expressed on activated CD4<sup>+</sup> and CD8<sup>+</sup> T cells, and is constitutively expressed on regulatory T cells (T<sub>regs</sub>). While GITR appears to act as a conventional co-stimulatory receptor in effector CD4<sup>+</sup> and CD8<sup>+</sup> T cells, the role of GITR on T<sub>regs</sub> remains controversial.

Here we utilized a phosphoproteomics approach to identify and compare GITR-mediated intracellular signaling events in induced-T<sub>regs</sub> (iT<sub>regs</sub>), CD4<sup>+</sup>, and CD8<sup>+</sup> cytotoxic T cells. The GITR agonist monoclonal antibody DTA-1 was used to stimulate the T cell subsets and GITR agonism was validated via NFκB pathway activation and phospho-JNK induction. PTMScan® immobilized metal affinity chromatography (IMAC) beads were used to enrich Ser/Thr/Tyr-modified peptides, and PTMScan® Phospho-Tyrosine Rabbit mAb (P-Tyr-1000) was used to capture an additional pool of phospho-Tyr modified peptides (Cell Signaling Technology, Inc.).

## CONCLUSIONS

- The phosphoproteomics data revealed distinct protein phosphorylation events between DTA-1-treated iT<sub>regs</sub>, CD4<sup>+</sup> T<sub>eff</sub>, and CD8<sup>+</sup> T cells.
- DTA-1 induced phospho-p38 MAPK pathway activation in iT<sub>regs</sub>, but not in CD4<sup>+</sup> T cells or CD8<sup>+</sup> cytotoxic T cells.
- Induction of p38 MAPK signaling is functionally important as inhibition of p38 abrogated GITR-induced proliferation of iT<sub>regs</sub>.
- These data demonstrate that GITR agonism induces unique phospho-proteome profiles between T cell subsets, and this differential signaling leads to unique responses that will need to be considered when targeting GITR with immunotherapeutic strategies.



**Schematic of phosphoproteomics analysis of anti-GITR treated T cell subsets**  
 Cell extracts from GITR agonist DTA-1 or isotype control treated iT<sub>regs</sub>, CD4<sup>+</sup> T<sub>eff</sub>, and CD8<sup>+</sup> T cells were digested into peptides, purified over reversed phase columns, enriched with pY antibodies or IMAC technology, and run on a LC-MS/MS. The data set was searched/score filtered, and relative quant data were generated. pY and IMAC enrichments were performed on independent sample preps.

## Methods

**Figure 1:**

## Results

**Table 1:**

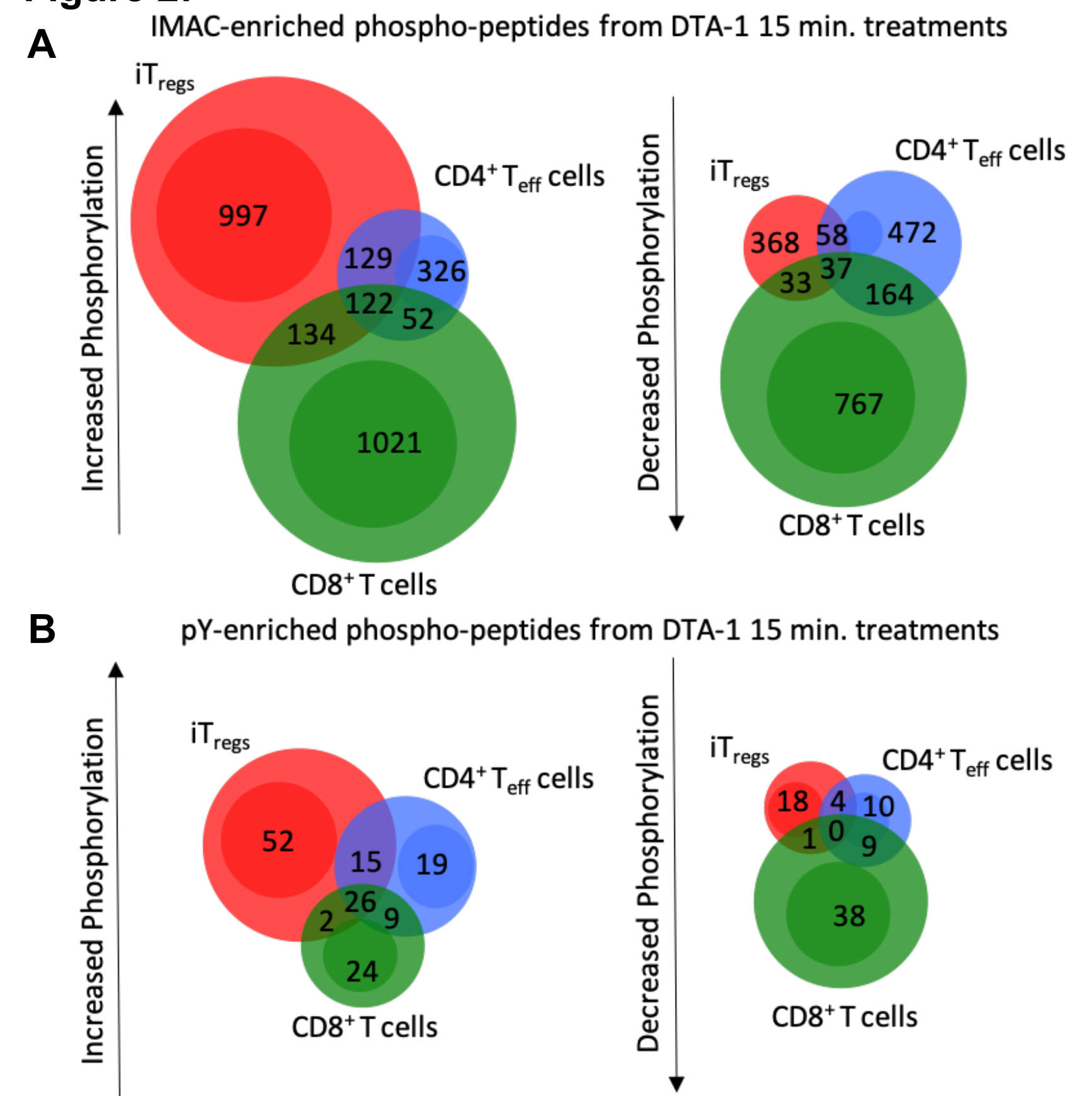
		IMAC	pY
Total Number of Detected Peptides		18,830	771
Total Number of Significant FC		5,645	244

iT <sub>regs</sub>		CD4 <sup>+</sup> T <sub>eff</sub> cells		CD8 <sup>+</sup> T cells	
Time Point	15 min.	Time Point	15 min.	Time Point	15 min.
Enrichment	IMAC	pY	Enrichment	IMAC	pY
Increase	1382	95	Increase	629	69
Decrease	496	23	Decrease	731	23
Increase			Increase	1,329	61
Decrease			Decrease	1,001	48

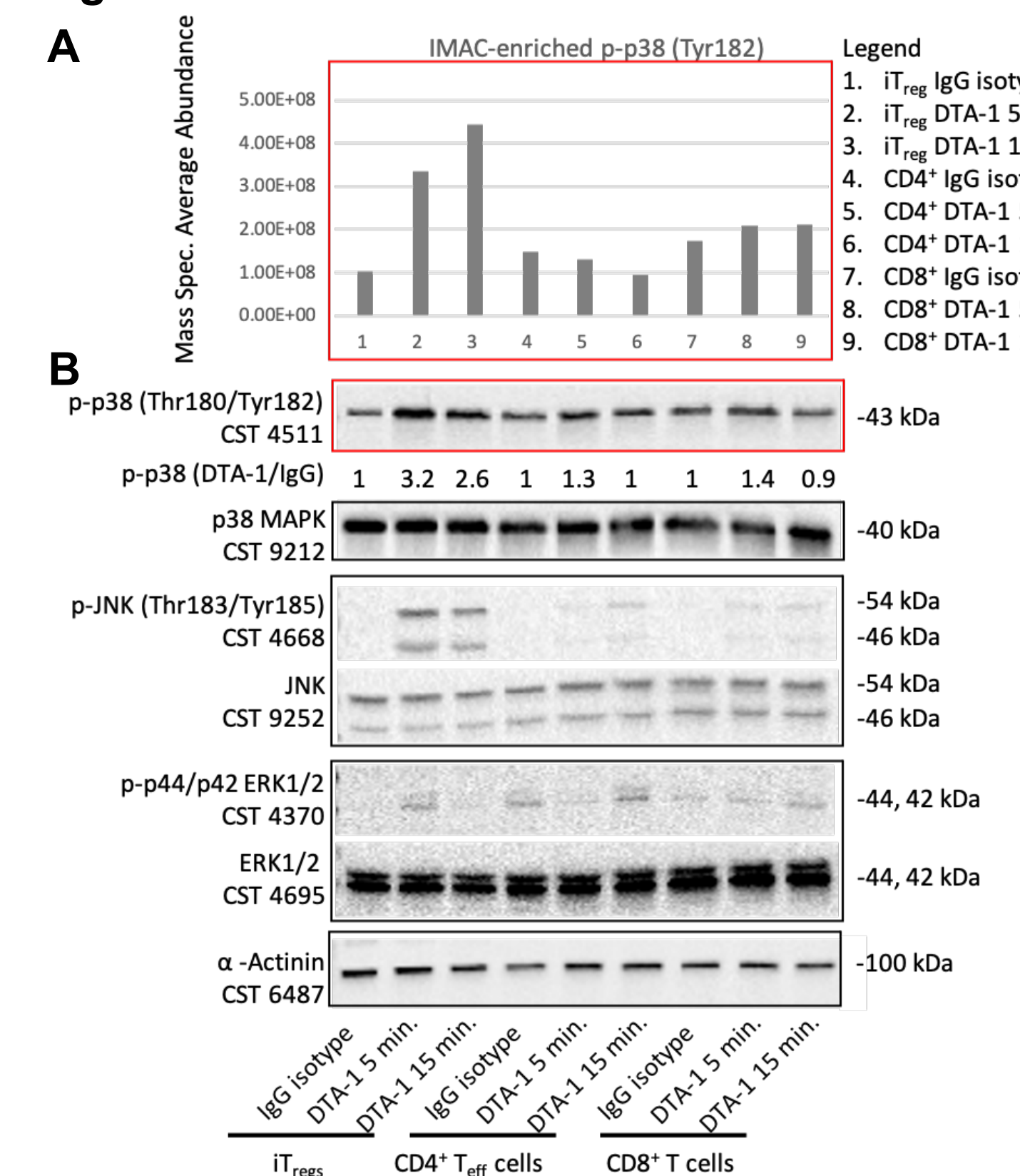
**Summary of anti-GITR treatment IMAC and pY Enrichment Results.** IMAC and pY-enriched phospho-peptide data from T cell subsets treated with 5 µg/mL IgG isotype control for 15 min. or 5 µg/mL DTA-1 for 5 or 15 min. Data are from two replicate LC-MS/MS runs per sample and were filtered to reflect nonredundant phosphorylation sites. The number of significant increases and decreases are shown for each T cell subset at 15 min. DTA-1 timepoints relative to IgG isotype control. Significant fold change increases are defined as the number of upregulated phospho-peptides ≥ 2.5 fold change and the number of decreases are defined as the number of downregulated phospho-peptides ≤ -2.5 fold change.

**Figure 2:**



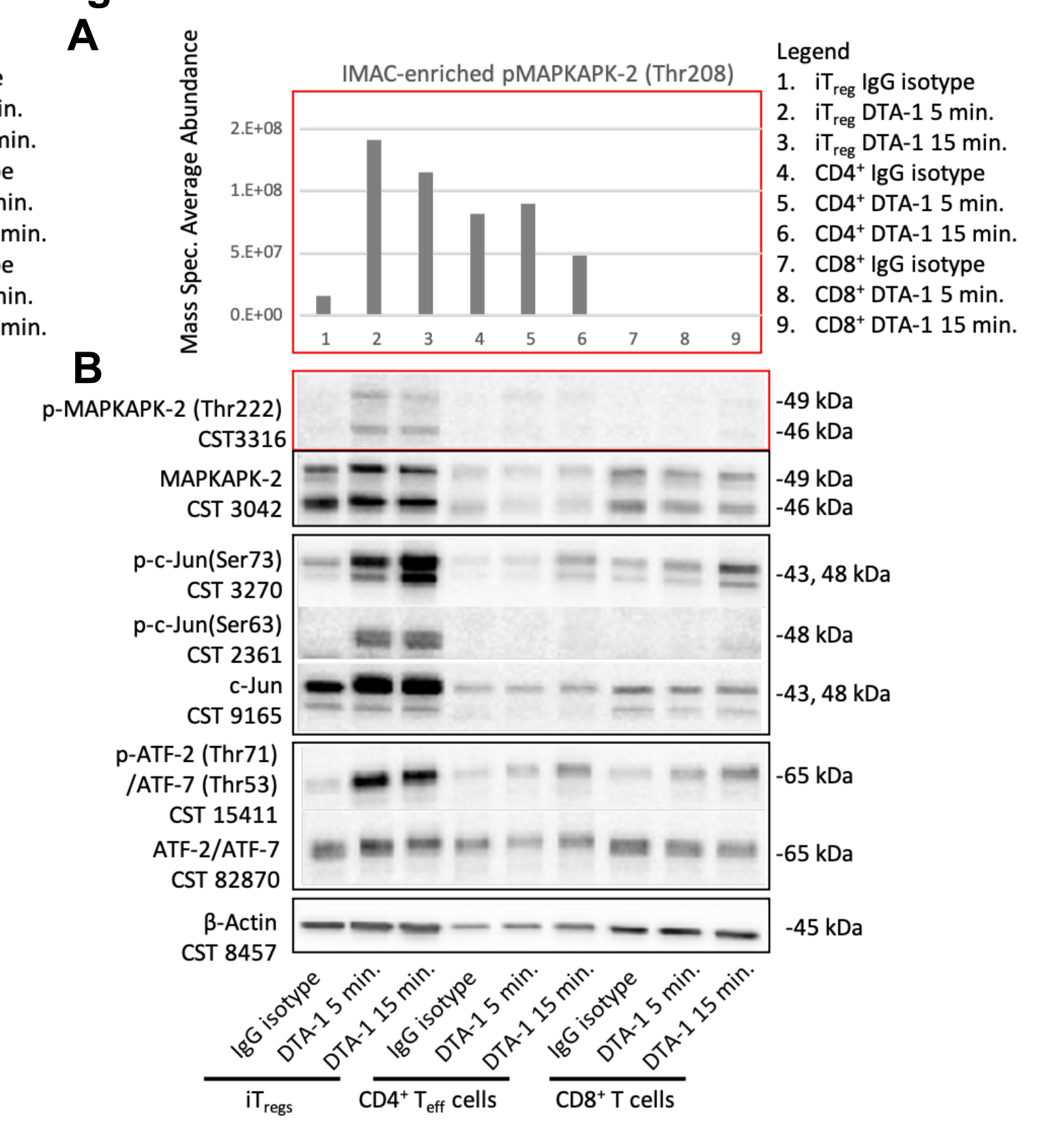
**Venn diagrams of DTA-1-regulated phosphorylation events comparing iT<sub>regs</sub>, CD4<sup>+</sup> T<sub>eff</sub>, and CD8<sup>+</sup> T cells.** (A) The number of unique IMAC-purified phospho-peptides that had elevated phosphorylation ≥ 2.5 fold (left), and reduced phosphorylation ≤ -2.5 fold change (right). (B) Venn diagrams showing the number of pY antibody-purified phospho-peptides that had elevated phosphorylation ≥ 2.5 fold change (left), and reduced phosphorylation ≤ -2.5 fold change (right). The relative size of the T cell subsets in the diagrams corresponds with the number of unique phospho-peptides detected. (A) is scaled for IMAC-purified peptides and can be compared across increases and decreases, and (B) is scaled for pY-purified peptides and can be compared across increases and decreases.

**Figure 3:**



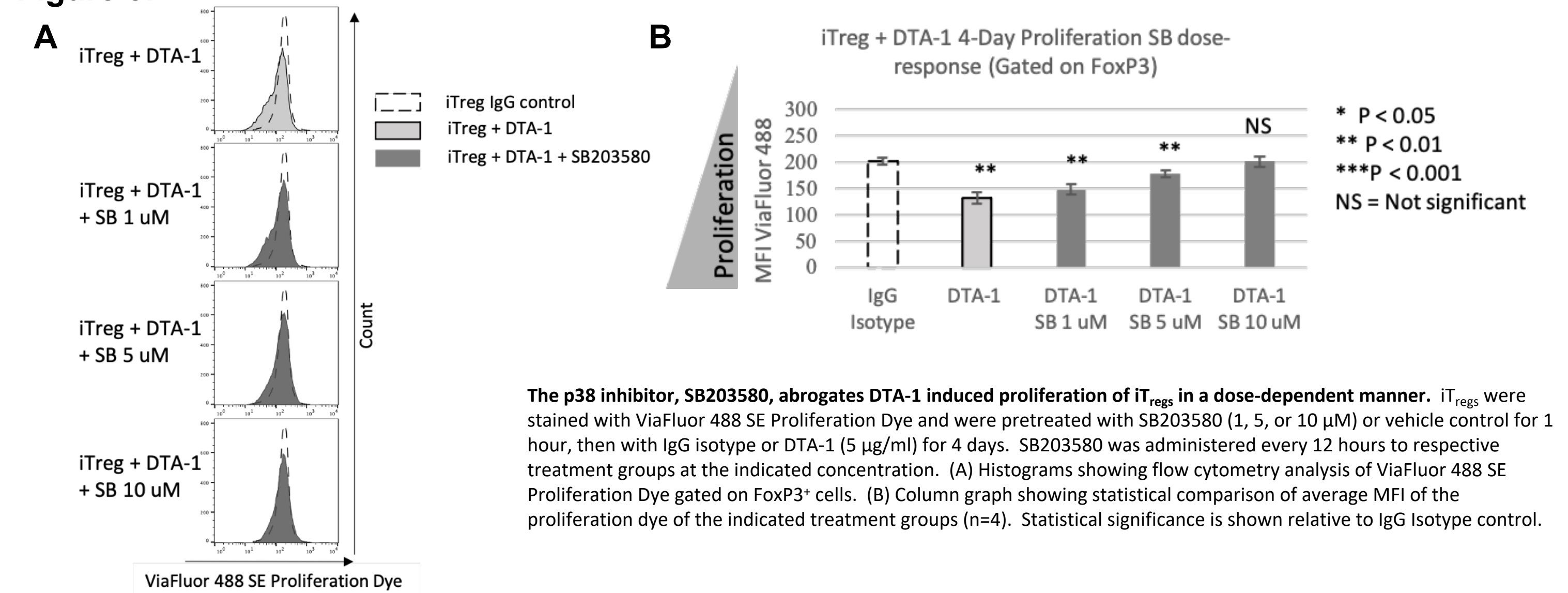
**GITR triggering induces p38 phosphorylation in iT<sub>regs</sub> but not in CD4<sup>+</sup> T<sub>eff</sub> cells and CD8<sup>+</sup> T cells.** (A) Mass spectrometry measured average abundance (intensity) of p-p38(Tyr182) detected in the T cell subsets. (B) Western blot analysis of the three main phospho and total MAPK signaling pathway members. P-p38 is highlighted in red boxes to easily compare phospho-proteomics and western blotting detection. Column layout of samples in (A) corresponds with western blot lanes in (B). Relative fold change of p-p38 is shown normalized to IgG control for each subset.

**Figure 4:**



**GITR triggering induces signaling downstream of p38 phosphorylation in iT<sub>regs</sub> differentially compared to CD4<sup>+</sup> T<sub>eff</sub> cells and CD8<sup>+</sup> T cells.** (A) Mass spectrometry measured average abundance (intensity) of p-MAPKAPK-2(Thr208) detected in the T cell subsets. P-MAPKAPK-2 (Thr222) is the human phospho site that the antibody product is named for, and (Thr208) is the mouse site. (B) Western blot analysis of phospho and total proteins downstream of p38.

**Figure 5:**



**The p38 inhibitor, SB203580, abrogates DTA-1 induced proliferation of iT<sub>regs</sub> in a dose-dependent manner.** iT<sub>regs</sub> were stained with ViaFluor 488 SE Proliferation Dye and were pretreated with SB203580 (1, 5, or 10 µM) or vehicle control for 1 hour, then with IgG isotype or DTA-1 (5 µg/ml) for 4 days. SB203580 was administered every 12 hours to respective treatment groups at the indicated concentration. (A) Histograms showing flow cytometry analysis of ViaFluor 488 SE Proliferation Dye gated on FoxP3<sup>+</sup> cells. (B) Column graph showing statistical comparison of average MFI of the proliferation dye of the indicated treatment groups (n=4). Statistical significance is shown relative to IgG Isotype control.