

Immunomodulation of Primary Human T Cell Activation Using High-throughput Screen of FDA-approved Small Molecule Drugs

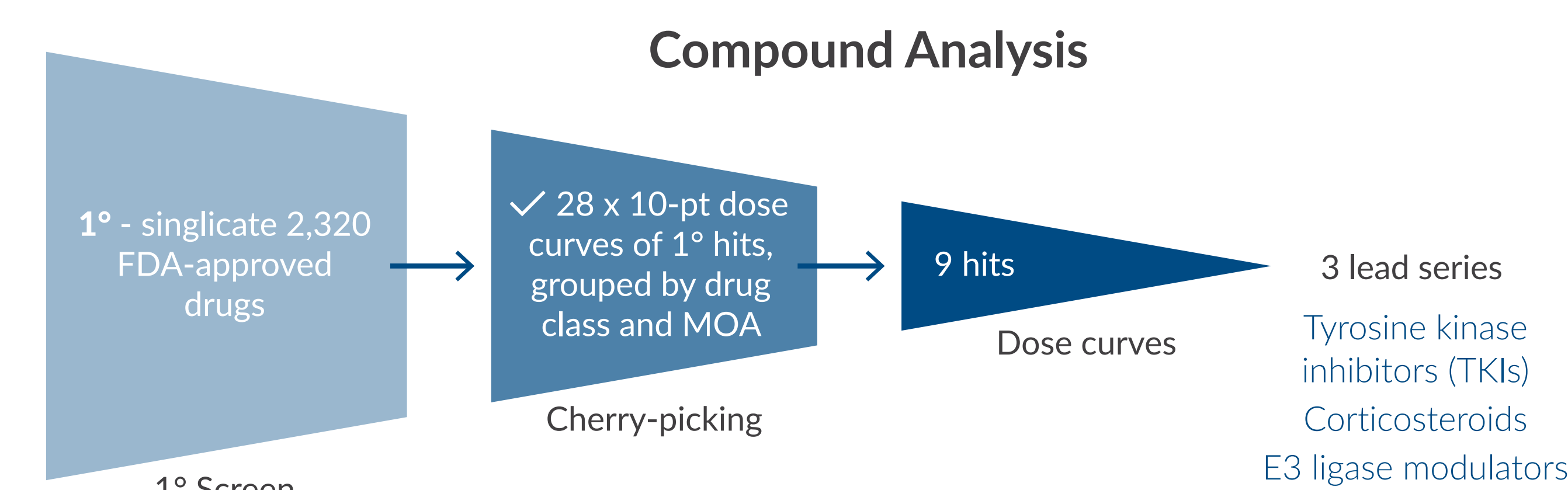
Marie Foss and David Hoffman
Cayman Chemical Company, Ann Arbor, MI

KEY FINDING By combining IL-2 secretion and viability assays, our screening platform accurately identifies multiple members of active compound classes and differentiates immunomodulatory or toxic activity for T cells.

ABSTRACT

Immunotoxicity and immunomodulation are frequent on- and off-target properties of therapeutics. The incorporation of immune function assays into structure-activity relationship (SAR) studies can provide valuable guidance for drug development, whether for drug safety and/or efficacy. In this high-throughput screen (HTS), we focus on the health and function of primary human T cells through utilization of an AlphaLISA IL-2 assay for IL-2 secretion and CellTiter-Glo® assay for cell survival and proliferation. When stimulated via the T cell receptor, T cells will begin expansion and secretion of the cytokine IL-2.

Incorporation of liquid handling automation allows for miniaturization of this primary cell-based assay to the 384-well format with improved throughput and reduction in material costs. Up to 28 × 10-point dose curves can be performed per 384-well plate in singlicate with replicates produced across multiple plates or 9 × 10-point dose curves performed in triplicate on a single 384-well plate. By utilizing cryopreserved peripheral blood mononuclear cells (PBMCs), assays can be regulated for batch-to-batch consistency and data turn around reduced to 48 hours from drug treatment.



HIGH-THROUGHPUT 1° SCREENING METHODS AND RESULTS

Our objective was to optimize our workflow for performance of a HTS of 2,320 FDA-approved small molecule drugs using an IL-2 AlphaLISA assay to detect changes in T cell activation. We sought to test whether modulation of IL-2 secretion by known drugs could be measured and characterized for future SAR studies. We developed workflows for both 1° screening at a single drug concentration, to be followed by cherry-picking of hits for dose curves and cytotoxicity evaluation, for identification of compounds that are cytotoxic versus immunomodulatory.

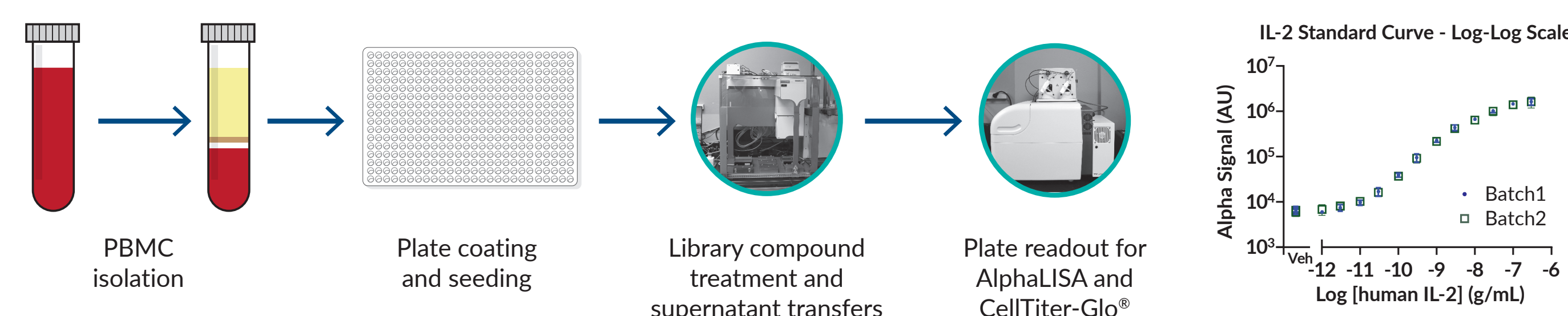


FIGURE 1 – Workflow for cell isolation, treatment, and readout.

- Isolate human PBMCs from whole blood.
- Activate T cells by seeding on αCD3/αCD28-coated 384-well plates.
- Treat with 2,320 FDA-approved compounds at 250 nM from 8 × 384-well library stock plates in DMSO.
- Staurosporine (SP) was a control for IL-2 inhibition and cytotoxicity.
- Incubate for 48 hours and transfer supernatant for IL-2 detection by AlphaLISA.

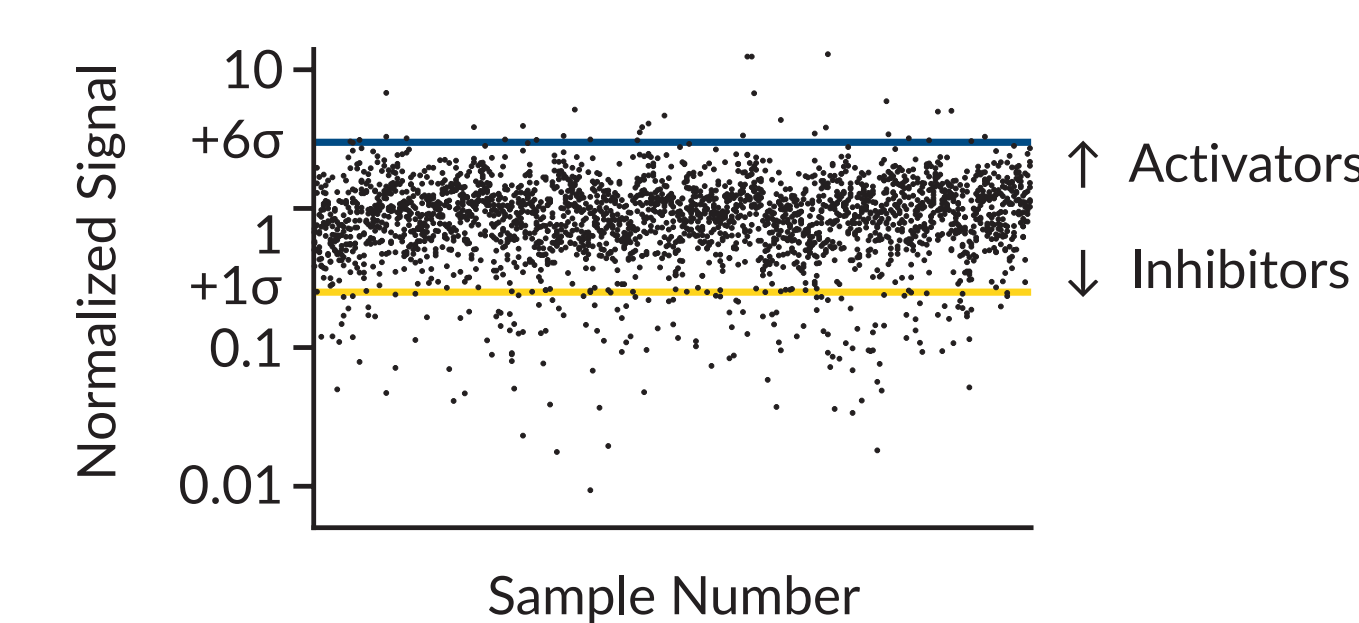


FIGURE 2 – Dot plot of 1° screen hits.

Cutoffs for IL-2 secretion activators were set relative to the vehicle control (0.025% DMSO) and the cutoff for inhibitors set relative to the negative control (250 nM SP). The sample values are normalized to the average signal across the full dataset.

Compound Class	Identified Members from 1° Screen Hits
Tubulin inhibitors	colchicine, vinorelbine, vinblastine
E3 ligase modulators	lenalidomide, pomalidomide
TKIs	dasatinib, ponatinib, masitinib, erlotinib, regorafenib, bosutinib, dovitinib, foretinib, ibrutinib, (R)-crizotinib (but not imatinib, nilotinib)
Corticosteroids	halobetasol, triamcinolone, fluocinonide, clobetasol, mometasone, dexamethasone
Anthracyclines	daunorubicin, idarubicin (but not epirubicin)

TABLE 1 – Drug classes identified by multiple hits in the 1° screen.

The HTS highlighted several classes of compounds that affected IL-2 secretion by T cells. Cutoff values for activation were set relative to DMSO vehicle controls and identified 1.0% of the FDA-approved library as IL-2 secretion stimulators at 250 nM treatment. Cutoff values for inhibition were set based on the negative control and identified 4.4% of the library as inhibitors of IL-2 production. IL-2 secretion activators are highlighted in blue and inhibitors are highlighted in yellow.

HIT PICKING FOR REPRODUCIBILITY AND EVALUATION OF POTENCY

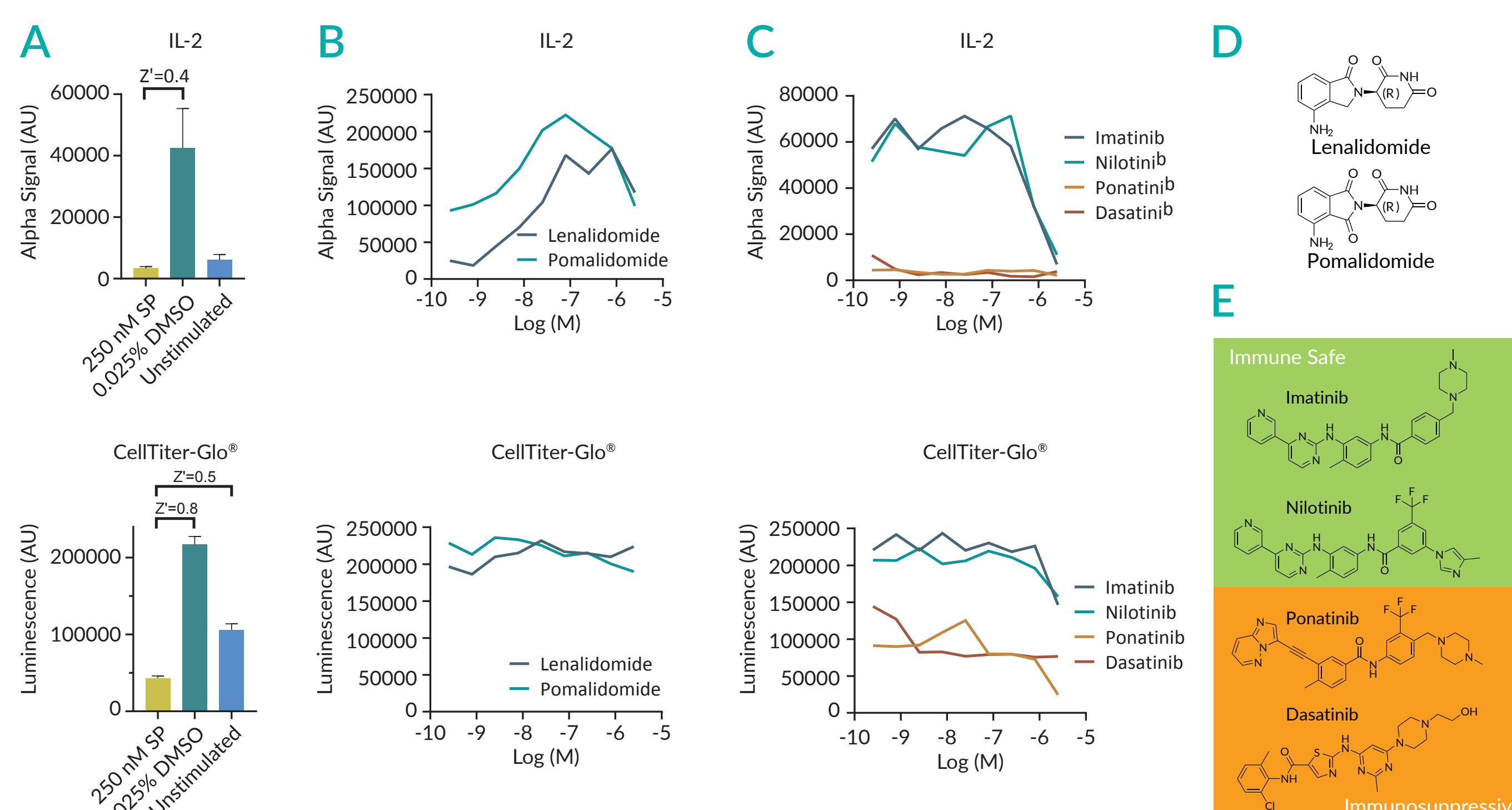


FIGURE 3 – Singlicate dose curves of cherry-picked hits based on drug class.

We verified the activity of 28 of the top activators and inhibitors grouped by drug class from the 1° screen to eliminate false positives. **A)** Control values from singlicate dose curve readouts. AlphaLISA Z' ranged from 0.4-0.6 during optimization and testing. CellTiter-Glo® Z' ranged from 0.7-0.8. **B)** From the E3 ligase modulators, the doses tested did not result in cell death. The effect at 250 nM in the 1° screen is reproducible. **C)** TKIs showed member-specific suppression of T cell activation, presenting opportunity for quantitative evaluation of SAR. **D)** E3 ligase modulator compound structures. **E)** TKI compound structures.

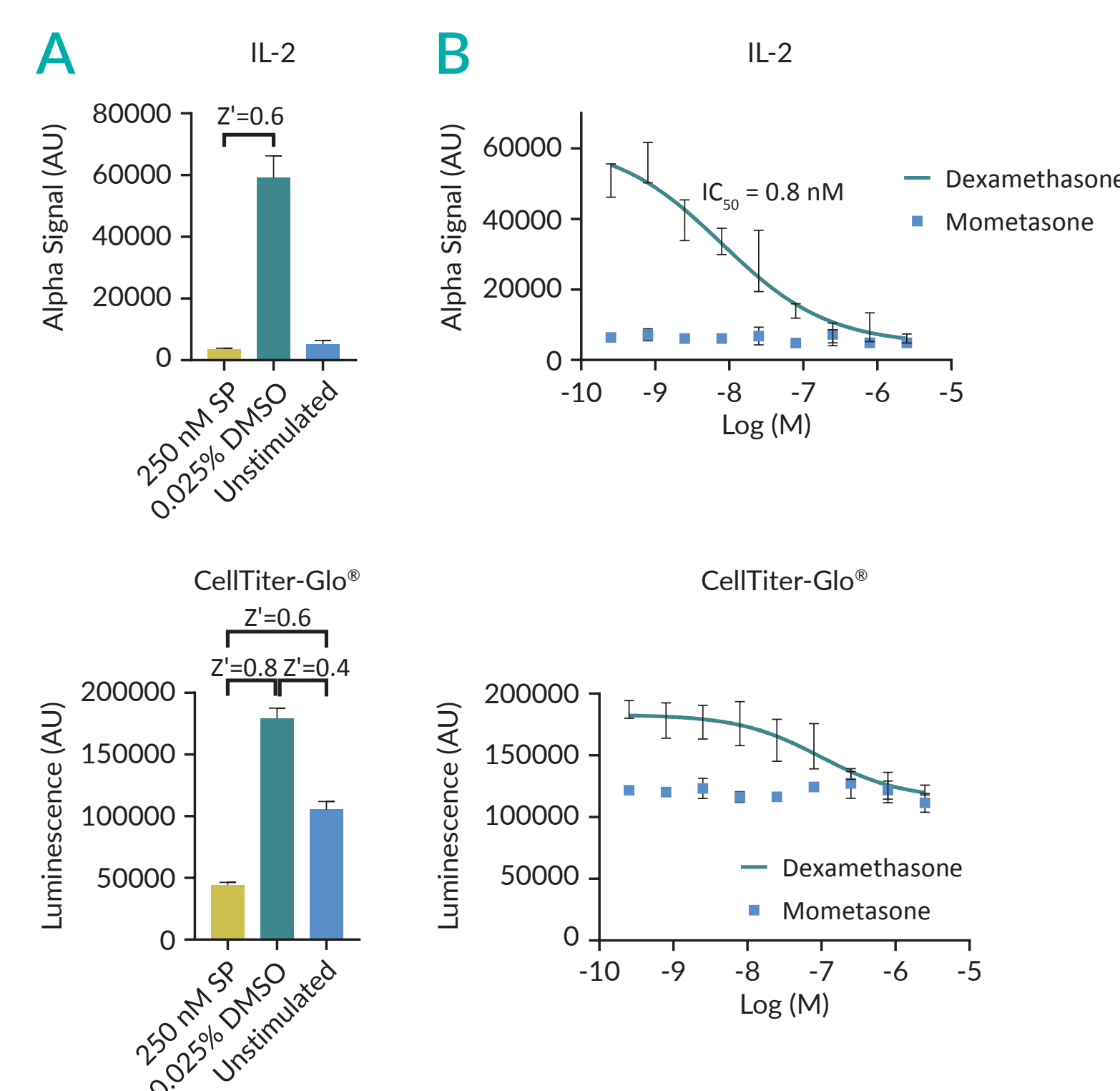


FIGURE 4 – Dose curves of verified hit compounds.

We performed dose curves in triplicate for nine compounds with verified activity in the 1° screen. **A)** Control values from singlicate dose curve readouts. **B)** We measured the inhibitory concentration at 50% inhibition (IC₅₀) for dexamethasone against IL-2 secretion at 0.8 nM. Typical values of glucocorticoid receptor activity for dexamethasone are low nM. Mometasone is known to bind the glucocorticoid receptor 22 times stronger than dexamethasone. **C)** Corticosteroid compound structures.

CONCLUSION

- The use of primary T cells gives a more direct *in vivo* comparison of small molecule drug activity for validation of mechanism of action for therapy.
- IL-2 secretion (T cell function) paired with cytotoxicity assays (T cell survival) can be used to differentiate small molecule immunotoxicity or immunomodulation of T cells, as well as direct cell killing mechanisms.
- T cell activators included certain tubulin inhibitors as well as E3 ligase modulators.²
- IL-2 suppressors that were not cytotoxic included corticosteroids and certain TKIs.
- Some TKIs have off-target effects on T cell receptor signaling at the tested dose, which highlighted immune safe members of this compound group that were tested during cherry picking for verification.³

References

- Mometasone furoate DrugBank DB14512. go.drugbank.com/drugs/DB14512.
- Gandhi, A.K., Kang, J., Havens, C.G., et al. Immunomodulatory agents lenalidomide and pomalidomide co-stimulate T cells by inducing degradation of T cell repressors Ikaros and Aiolos via modulation of the E3 ubiquitin ligase complex CRL4^{CRBN}. *BJ Haem.* **164**(6), 811-821 (2014).
- Lamble, A.J., Kosaka, Y., Laderas, T., et al. Reversible suppression of T cell function in the bone marrow microenvironment of acute myeloid leukemia. *Proc. Natl. Acad. Sci. USA.* **117**(25), 14331-14341 (2020).



Download this poster & learn more about our high-throughput screening services