

ATCC® 100 YEARS

Abstract #3959

Abstract

Background

[illegible]

Results

A Luminescence intensity HCC827 GAS-Luc2

IFN- γ concentration (ng/mL)	Fold Change
0	~10
0.01	~20
0.1	~30
1	~100
10	~180
100	~240
1000	~250

B Luminescence intensity HCC827 GAS-Luc2

Condition	Fold Change
Non-activated	~60
Activated	~100

C Luminescence intensity MG-63 GAS-Luc2

IFN- γ concentration (ng/mL)	Fold Change
0	~10
0.01	~10
0.1	~10
1	~20
10	~90
100	~170
1000	~230

D Luminescence intensity MG-63 GAS-Luc2

Condition	Fold Change
Non-activated	~5
Activated	~35

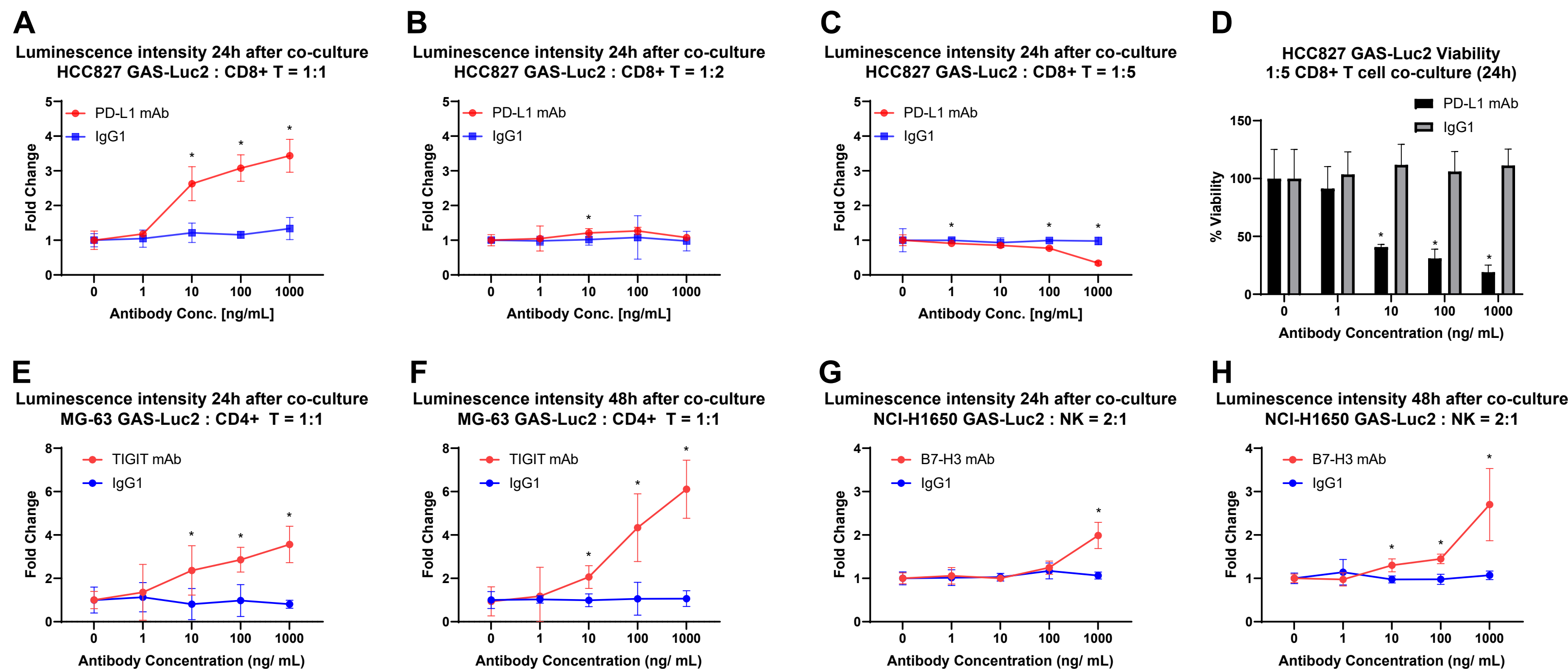
E Luminescence intensity NCI-H1650 GAS-Luc2

IFN- γ concentration (ng/mL)	Fold Change
0	~10
0.01	~10
0.1	~10
1	~30
10	~80
100	~130
1000	~140

F Luminescence intensity NCI-H1650 GAS-Luc2

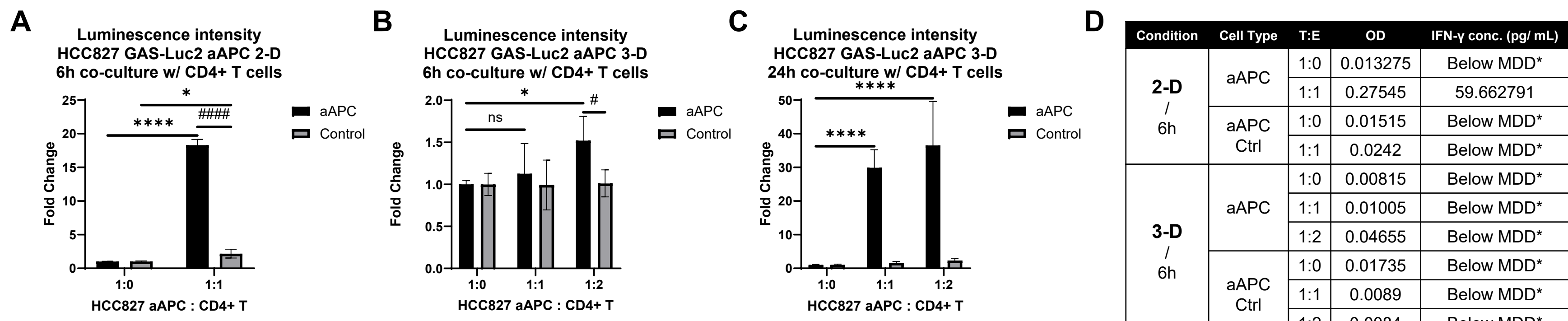
Condition	Fold Change
Non-activated	~5
Activated	~50

Co-culture with adaptive or innate immune cells yields luciferase expression in GAS-Luc2 cells

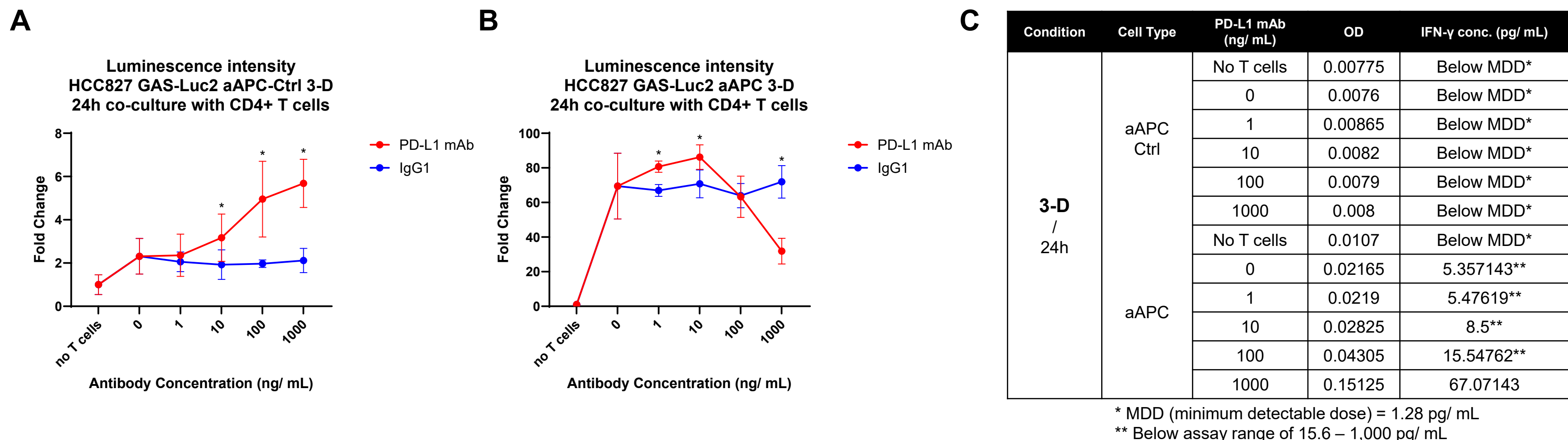


Key References

- ### **GAS-Luc2 cells demonstrate superior sensitivity to ELISA in 2-D and 3-D co-culture with T cells**



Comparison of GAS-Luc2 cells and ELISA in immune checkpoint inhibitor evaluation



Conclusion

- ATCC**

10801 University Boulevard, Manassas, Virginia 20110-2209

Phone: 800.638.6597

Email: sales@atcc.org

Web: www.atcc.org

© 2025 American Type Culture Collection. ATCC product identifiers marked with the TM symbol are trademarks owned by the American Type Culture Collection. Bright-Glo and Promega are trademarks or registered trademarks of Promega Corporation. SpectraMax and Molecular Devices are registered trademarks of Molecular Devices, LLC. Quantikine is a registered trademark of Research & Diagnostic Systems, Inc. R&D Systems is a registered trademark of Bio-Techne Corporation.