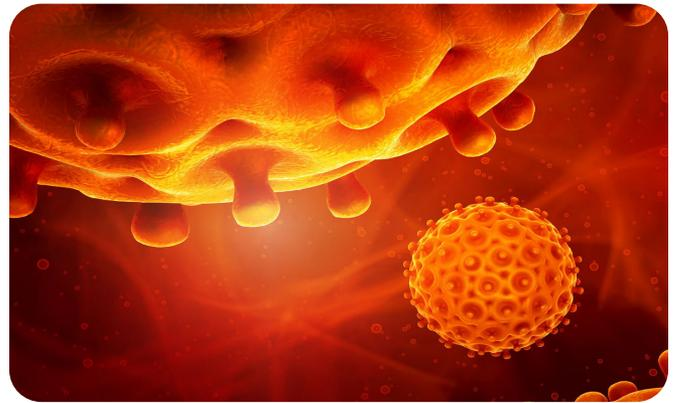
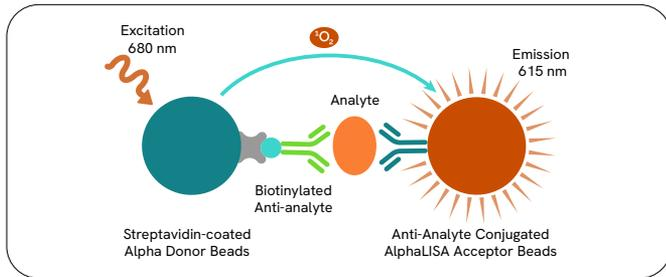




Innovative solutions in virology research.

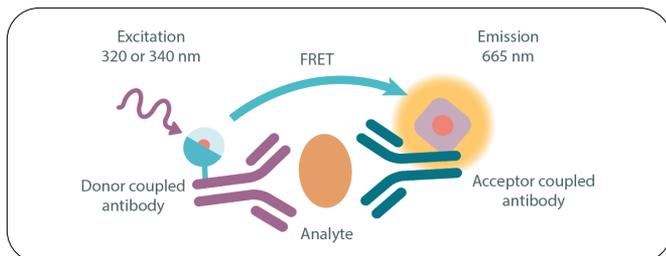
Alpha™ technology

Alpha is a versatile, bead-based platform that enables you to assay the most complex samples in one well and with no wash steps.



HTRF assays

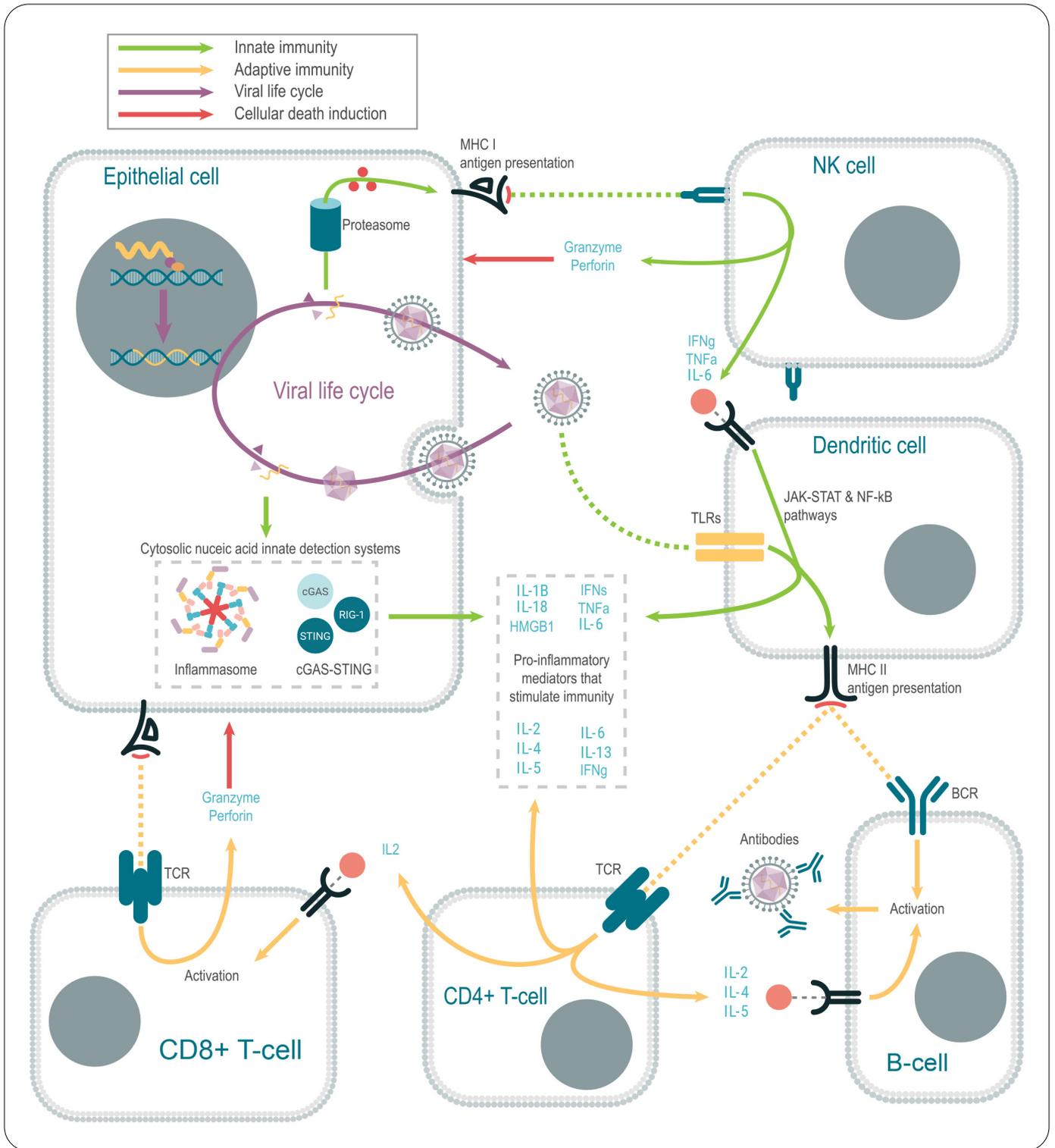
Our HTRF™ assays are fast, sensitive, homogeneous, and ready to use with no wash steps.



Revvity also provides luminescence assays based on firefly luciferase.

- Britelite™ plus, neolite™ and steadylite plus high to ultra high sensitivity reporter gene assay systems
- Cytotoxicity/cell proliferation assays based on ATP monitoring to assess the cytotoxic, cytostatic and proliferative effects of a wide range of drugs, biological response modifiers and biological compounds (ATPlite technology).

Antiviral immune response and immunization strategies



Available products

Table 1. Viral detection assay.

Virus quantification AlphaLISA Kits 500 test	Part #
HIV p24 (high sensitivity)	AL291C
Hepatitis B Virus e Antigen (HBeAg)	AL308HV
Hepatitis B Virus s Antigen (HBsAg)	AL3083HV
HCV cAg (hepatitis C virus core antigen)	AL3092HV
Influenza A Nucleoprotein	AL3084HV
Viral double-stranded RNA detection products	Part #
Viral double-stranded RNA detection Kit 500 test	64RNAPEG
Viral double-stranded RNA detection Kit standard	64RNATDA
Lysis buffer 16 mL	64B4FDD
Virus detection HTRF Kits 500 tests	Part #
SARS-CoV-2 nucleocapsid	63ADK113PEG
SARS-CoV-2 Spike S1 protein	63ADK114PEG
SARS-CoV-2 S1/ACE2 binding	64BDS1PEG
Virus detection AlphaLISA Kits, 500 tests	Part #
SARS-CoV2 Spike S1	AL3142C
SARS-CoV2 Nucleocapsid	AL3141C
SARS-CoV2 ACE2/S1	AL3146C

Table 2. Key cytokines.

Key Cytokines 500 tests	HTRF	AlphaLISA
IL-1b	62HIL1BPEG	AL3160C
IL-2	62HIL02PEG	AL3155C*
IL-6	62HIL06PEG	AL3025C
IL10	62HIL10PEG	AL3159C
IL-13	62HIL13PEG	AL240C
IL-18	62HIL18PEG	AL559C
HMGB1	62HMGPEG	
TNF- α	62HTNFAPEG	AL3157C*
IFN- α	62HIFNAPEG	AL576C
IFN- β	62HIFNBPEG	AL265C
IFN γ	62HIFNGPEG	AL3153C*
IFN- λ 2		AL3031C
TGF β -1	62HTGFBPEG	AL361C

*Available in biotin-free format

Table 3. COVID-19 cytokine storm.

COVID-19 cytokine storm assays 500 tests	HTRF	AlphaLISA
IL-6	62HIL06PEG	AL3025C
C Reactive Protein		AL233C

Table 4. Innate antiviral immunity.

Pathways	Assay 500 tests	HTRF	AlphaLISA SureFire Ultra or AlphaScreen SureFire
TLR	IRAK3	63ADK101PEG	
	IRAK4	63ADK108PEG	AL3117C
	JNK Phospho- T183/S185*	64JNKPEG	TGRJS500
	c-Jun Phospho- S63		TGRCJS500
	c-Jun Phospho- S73		TGRCJ2S500
	p38 MAPK Phospho- T180/Y182*	64P38PEG	ALSU-PP38-B500
cGAS-STING	Mouse STING phospho S365	64MSTGPEG*	
	cGAS total	64CGASTPEG	
	STING Phospho- S366*	64P62S4PEG	
	Human STING WT binding Kit	64BDSTGPEG	AL3144C
	Mouse STING WT binding Kit	64BDSTGMPEG	
	Human AQ STING binding Kit	64BDSTGQPEG	
	Human H232 STING binding Kit	64BDSTGHPEG	
	p62		AL3097C
TLR/ cGAS-STING/ Inflammasome/ NF-kB	TBK1 Phospho- S172*	64TBKPEG	
	IRF3 Phospho- S385/386	6FRF3PEG	ALSU-PIRF3-A500*
	IKBA phospho S32/36		ALSU-PIKBA-A500*
	Ikkα Phospho- S176/180	64KKAS6PEG*	ALSU-PIKKA-A500
	Ikkβ Phospho- S177/181	64KKBPEG	ALSU-PIKKB-A500*
	NF-kB	64NFBPEG	ALSU-PNFKB-A500
JAK-STAT	Human Phospho-IkBα (Ser32/36) Detection Kit*		ALSU-PIKBA-A500
	Mouse Phospho-IkBα (Ser32/36) Detection Kit*		ALSU-PIKBA-A500
	Human Phospho-IKKβ (Ser177/181) Detection Kit*		ALSU-PIKBA-A500
	RIG-I Binding Kit	64BDRIGIPEG	
	STAT1 Phospho- Y701*	63ADK026PEG	ALSU-PST1-A500
	STAT1 Phospho- S727*		ALSU-PST1-B500
	STAT3 Phospho- Y705*	62AT3PEG	ALSU-PST3-A500
	STAT4 Phospho- Y693		ALSU-PST4-A500
	STAT5 Phospho- Y694	64AT5PEG	
	STAT5 Phospho- Y694/699*		ALSU-PST5-B500
	STAT6 phospho Y641		ALSU-PST6-A500
	JAK1 Phospho-Tyr1034/1035 Detection Kit*	64JAK1Y10PEG	
	JAK2 Phospho-Tyr1007/1008 Detection Kit*	64JAK2Y10PEG	
JAK3 Total Detection Kit*	64JAK3TPEG		

*Available in Phospho-- and total protein format

Table 5. Viral neutralization assays.

Reporter gene reagents	PART #
Britelite Plus*	606676*
neolite*	601671*
steadylite Plus*	606675*

*Exists in 10/100/1,000 mL

Table 6. Cytotoxicity and viability assays.

Luminescence assays	PART #
ATPlite 1step	6016736*
ATPlite	6016943**

*Exists in 10/100/1,000 mL; **Exists in 300/1,000/5,000/10,000 tests

Table 7. Viral neutralization assays.

Fluorescence assays 500 tests	HTRF	Alpha
pH2AX	64H2XPEG	ALSU-PH2X-A500
Cleaved PARP	64PARPEG	

Table 8 (A). Adaptive antiviral immunity.

T-cell activation pathway 500 tests	HTRF	AlphaLISA SureFire Ultra or AlphaScreen SureFire
CD3 ζ Phospho-Tyr142 Detection Kit*		ALSU-PCD3-B500*
CD3 ζ Phospho-Tyr83 Detection Kit		ALSU-PCD3-A500*
ZAP70 Phospho- Y319*		ALSU-PZAP-A500
SLP76 Phospho- S376*	63ADK076PEG	ALSU-PSLP-A500
Human SLP76 Phospho-S376/total SLP76 Detection Kit		MPSU-PTSLP-K500
Mouse SLP76 Phospho-S376/total SLP76 Detection Kit		MPSU-PTSLP-K500
SLP76 phospho Y145 (mouse)		ALSU-PSLP-E500
SLP-76 Phospho-T145 Detection Kit	64SLPY5PEG	
ERK1/2 Phospho- T202/Y204*	64AERPEG	ALSU-PERK-A500

Table 8 (A). Adaptive antiviral immunity (Contd.)

T-cell activation pathway 500 tests	HTRF	AlphaLISA SureFire Ultra or AlphaScreen SureFire
CREB Phospho- S133*	64CREPEG	ALSU-PCREB-A500
SHP1 Phospho- Y536*		ALSU-PSHP1-A500
SHP1 Phospho- Y564*	64SH1PEG	ALSU-PSHP1-B500
SHP2 Phospho- Y542*	64SH2PEG	ALSU-PSHP2-A500
SHP2 Phospho- Y580*		ALSU-PSHP2-B500
JNK Phospho- T183/S185*	64JNKPEG	
c-Jun Phospho- S63		TGRCJS500
c-Jun Phospho- S73		TGRCJ2S500
NF-kB p65 Phospho- S536*	64NFBPEG	ALSU-PNFKB-A500
mTOR Phospho- S2448	64TORPEG	ALSU-PMTOR-C500
mTOR Phospho- S2481		ALSU-PMTOR-B500

*Available in Phospho-- and total protein format

Table 8 (B). Adaptive antiviral immunity.

B-cell activation pathway 500 tests	HTRF	AlphaLISA SureFire Ultra or AlphaScreen SureFire
SYK Phospho- Y352*		ALSU-PSYK-B500
SYK Phospho- Y525/526*	64SYKY525PEG	ALSU-PSYK-A500*
BTK Phospho- Y223*	63ADK017PEG	ALSU-PBTK-A500
BTK Phospho- Y551*		ALSU-PBTK-B500
ERK1/2 Phospho- T202/Y204*	64ERKPEG	ALSU-PERK-A500
JNK Phospho- T183/S185*	64JNKPEG	TGRJS500
c-Jun Phospho- S73		TGRCJ2S500
NF-kB p65 Phospho- S536*	64NFBPEG	ALSU-PNFKB-A500

*Available in Phospho-- and total protein format

Table 9. Selected KITS and reagents for immunoglobulin detection.

AlphaLISA detection KITS	AlphaLISA Acceptor Beads	HTRF KITS	HTRF labeled antibodies
IgM (bovine)	Anti-Bovine IgA Acceptor Beads	Human Fc	PAb Anti Human IgG-Eu cryptate
IgG1(bovine)	Anti-Bovine IgG Acceptor Beads	Human IgG	PAb Anti Human IgG-Tb cryptate
IgG2 (bovine)	Anti-Bovine IgG1 Acceptor Beads	Human IgG Kappa	Pab Anti Human IgG-d2
IgE (mouse)	Anti-Bovine IgG2 Acceptor Beads	Human IgG Lambda	PAb Anti Human IgG-XL665
IgG (human)	Anti-Bovine IgGM Acceptor Beads	Mouse Fc	MAB Anti Human Kappa-Eu cryptate
IgG (pan mouse)	Anti-Chicken IgY Acceptor Beads	Rabbit Fc	MAB Anti Human Kappa-d2
IgG Fc (human)	Anti-Goat (Fc-specific) Acceptor BeadsT		PAb Anti Mouse IgG-Eu cryptate
IgG (mouse)	Anti-Human (Fc-specific) Acceptor Beads		PAb Anti Mouse IgG-Tb cryptate
IgG1 (human)	Anti-Human IgE Acceptor Beads		Pab Anti Mouse d2
IgG1 (mouse)	Anti-Human (Fab specific) IgG Acceptor Beads		PAb Anti Mouse IgG-XL665
IgG2a (mouse)	Anti-Human IgG1 Acceptor Beads		PAb Anti Rabbit IgG-Eu cryptate
IgG2b (mouse)	Anti-Human IgG2 Acceptor Beads		PAb Anti Rabbit IgG-Tb cryptate
IgG3 (mouse)	Anti-Human IgG4 Acceptor Beads		PAb Anti Rabbit IgG-d2
IgG4 (human)	Anti-Human IgGE Acceptor Beads		PAb Anti Rabbit IgG-XL665
IgM (human)	Anti-Mouse (Fc-specific) IgG Acceptor Beads		Protein A-Eu cryptate
IgM (mouse)	Anti-Mouse IgG Acceptor Beads		Protein A-XL665
Kappa light chain	Anti-Mouse IgG1 Acceptor Beads		
	Anti-Mouse IgG2a Acceptor Beads		
	Anti-Mouse IgG2b Acceptor Beads		
	Anti-Mouse IgG3 Acceptor Beads		
	Anti-Mouse IgGM Acceptor Beads		

Custom labeling and custom assay development services

Don't see the assay you need listed here? Revvity has partnered with industry and academia professionals to develop custom solutions to dramatically shorten the drug development process using our powerful, accurate HTRF™,

DELFIATM and AlphaLISATM technologies. An experienced Project Manager will evaluate your project's chance of success, then work with you to create a customized assay for the target you need to advance your CAR-T research.

