



revvity

# SPA from Revvity: a clear advantage over other scintillation technology.

Combine the sensitivity of radioactive assays with the convenience of a homogeneous (no wash steps) assay for clearer research results using scintillation proximity assay (SPA) technology.

Revvity's SPA bead-based platform, consisting of imaging and scintillation beads, is a versatile, convenient, and cost-effective technology for use across various applications, including enzymatic activity assays, receptor-ligand binding studies, radioimmunoassays, and molecular interactions.

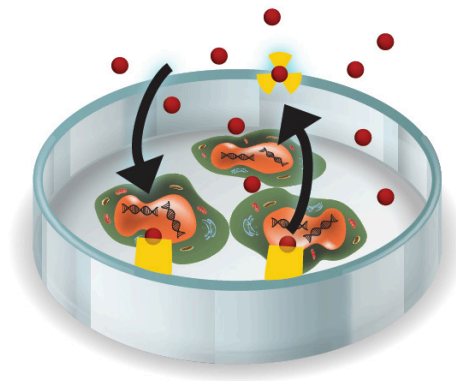


Scintillation-coated CytoStar-T™ plates are ideal for *in situ* cellular analysis, receptor binding assays, cell proliferation and viability, and drug metabolism studies. The advantages of SPA from Revvity have never been clearer.

CytoStar-T scintillating microplates: radiometric detection in whole cells in real time

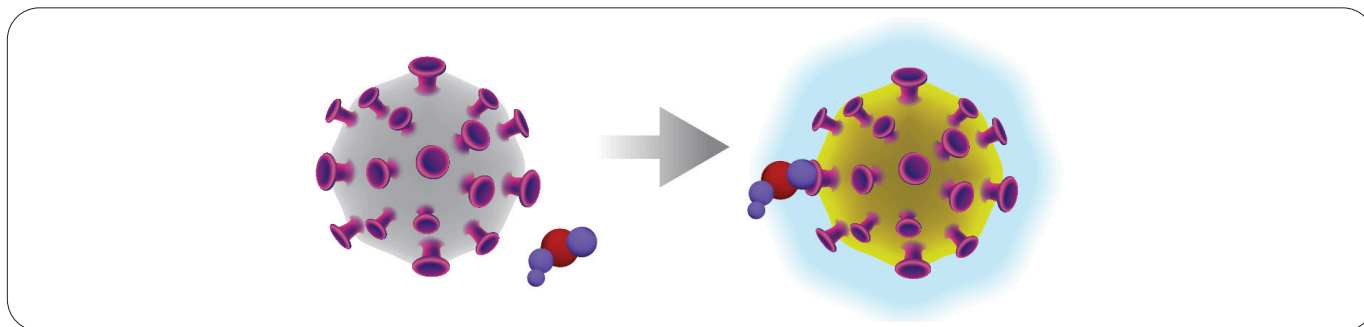


CytoStar-T 96 well plate



Microplate	Pack size	Product code
CytoStar-T - 96 well	5 pack	RPNQ0162
	100 pack	RPNQ0163
CytoStar-T - 384 well	5 pack	RPNQ0165
	100 pac	RPNQ0166

SPA scintillation beads: blue light emitting beads read on PMT-based radiometric detectors\*



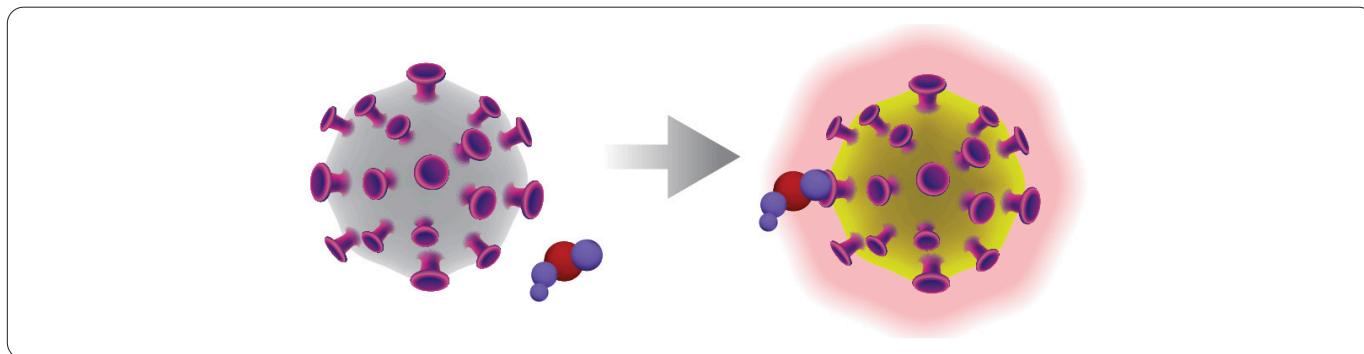
Coating	Application	Core bead type	Pack size	Product code
Streptavidin	Capture of biotinylated proteins or peptides for use in enzyme assays or molecular interaction studies	PVT	50 mg 150 mg 500 mg 2G 25 x 2G	RPNQ0006 RPNQ0009 RPNQ0007 RPNQ0066 RPNQ0067
		YSi	75 mg 250 mg	RPNQ0015 RPNQ0012
Wheat germ agglutinin (WGA)	Binds cell membranes and internal components for receptor binding studies	PVT	100 mg 500 mg 25 x 500 mg 2G	RPNQ0252 RPNQ0001 SPQ0031 RPNQ0060
		YSi	250 mg 1G	RPNQ0011 RPNQ0023
WGA polyethyleneimine (PEI) type A	Extra coating for reduced non-specific binding	PVT	500 mg 2G 25 x 2G	RPNQ0003 RPNQ0061 RPNQ0064
WGA PEI type B	Extra coating for reduced non-specific binding	PVT	500 mg 25 x 2G	RPNQ0004 RPNQ0065
PEI	Extra coating for reduced non-specific binding	PVT	500 mg	RPNQ0097
Poly - L- lysine	Enhances binding of negatively charged membranes	YSi	1000 mg	RPNQ0010
PDE	Measure of phosphodiesterase activity	YSi	500 mg 2G	RPNQ0150 RPNQ0024
RNA binding beads	YSi beads that interact with primary Phosphate groups in nucleotides, oligos, RNA and DNA; membranes can also bind	YSi	500 mg	RPNQ0013
Copper chelate	Capture and assay of His-tag fusion proteins or their binding partners	PVT	250 mg	RPNQ0095
		YSi	125 mg	RPNQ0096

Coating	Application	Core bead type	Pack size	Product code
Glutathione	Capture and assay of GST-fusion proteins	PVT	750 mg	RPNQ0030
		YSi	50 mg	RPNQ0033
Protein A	Binding of antibodies via the Fc portion of the antibody for RIA	PVT	500 mg	RPNQ0019
		YSi	500 mg 2G	RPN143 RPNQ0068
Anti-mouse antibody	Capture of secondary antibody	PVT	500 mg	RPNQ0017
		YSi	500 mg	RPN141
Anti-rabbit antibody	Capture of secondary antibody	PVT	500 mg	RPNQ0016
		YSi	500 mg	RPN140
SPA Scintillation Select-a-Bead Kit	Receptor binding studies	PVT & YSi	100 mg of 5 types**	RPNQ0250

\* Such as Revvity MicroBeta<sup>2</sup>™

\*\* Contains 100 mg each of WGA PVT; WGA YSi; WGA PEI type A; WGA PEI type B; and Poly-L-Lysine to allow quick and convenient receptor assay development using SPA Scintillation Beads.

### SPA imaging beads: red light emitting beads read on CCD-based radiometric detectors\*



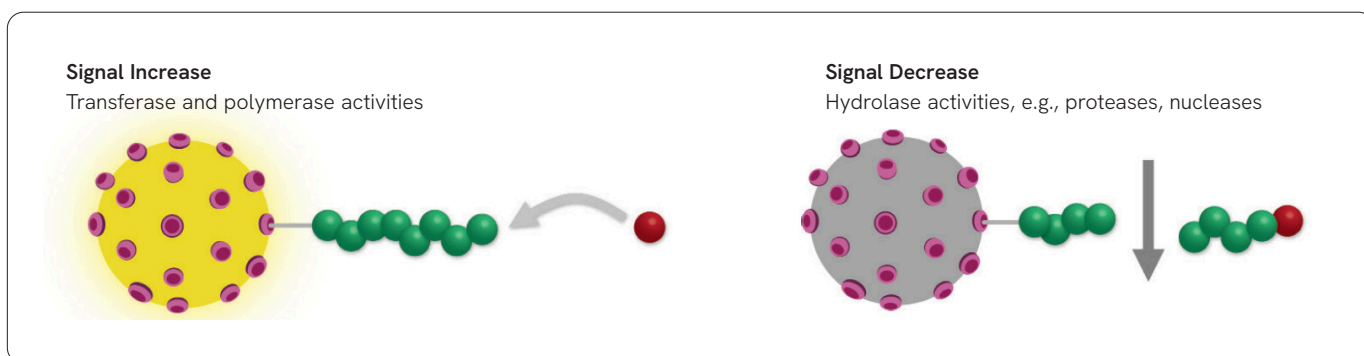
Coating	Application	Core bead type	Pack size	Product code
Streptavidin	capture of biotinylated proteins or peptides for use in enzyme assays or molecular interaction studies	PS	50 mg 500 mg	RPNQ0263 RPNQ0261
		YOx	50 mg 500 mg	RPNQ0096 RPNQ0271
Wheat germ agglutinin (WGA)	binds cell membranes and internal components for receptor binding studies with partially purified cell membrane preps or fractionated, solubilized receptor preps by immobilized receptors attached via glycosylation sites	PS	50 mg 500 mg	RPNQ0262 RPNQ0260
		YOx	50 mg	RPNQ0272
WGA polyethyleneimine (PEI) type A	PEI treatment decreases non-specific binding to beads	PS	50 mg 500 mg	RPNQ0286 RPNQ0287

Coating	Application	Core bead type	Pack size	Product code
PEI	PEI treatment decreases non-specific binding to beads	PS	50 mg 500 mg	RPNQ0297 RPNQ0098
		YOx	50 mg 500 mg	RPNQ0096 RPNQ0271
Poly - L- lysine	enhances binding of negatively charged membranes	PS	50 mg	
		YOx	2G	RPQ0328
Nickel chelate	capture and assay of His-tag fusion proteins	PS	500 mg	RPNQ0266
Glutathione	binding and assaying GST-fusion proteins	YOx	500 mg	RPNQ0278
Anti-mouse antibody	Affinity purified IgG designed for capture of secondary antibody	PS	500 mg	RPNQ0298
		YOx	500 mg	RPNQ0300
Anti-rabbit antibody	Affinity purified IgG designed for capture of secondary antibody	PS	500 mg	RPNQ0299
		YOx	500 mg	RPNQ0301
SPA Imaging Select-a-Bead Kit	Receptor binding studies	PS & YOx	500 mg of 6 types**	RPNQ0291

\* Such as the ViewLux™ CCD-based radiometric detector.

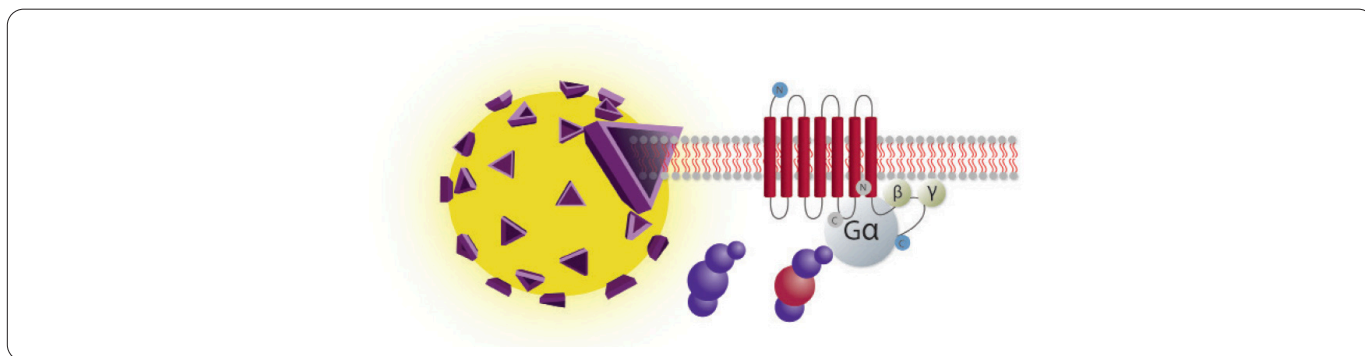
\*\* Contains 500 mg each of WGA PS; WGA YOx; WGA PEI type A; WGA PEI type B; PEI PS; and Poly-L-Lysine to allow quick and convenient receptor assay development using SPA Imaging Beads.

## SPA enzyme assays: radiometric detection of enzyme activity on bound target



Assay	Description	Pack size	Product code
PDE [ <sup>3</sup> H] cAMP assay	Phosphodiesterase assay	500 assays	TRKQ7090
PDE [ <sup>3</sup> H] cGMP assay	Phosphodiesterase assay	500 assays	TRKQ7100

## SPA receptor assays: radiometric detection of receptor-ligand binding



Assay	Description	Pack size	Product code
GTP $\gamma$ <sup>35</sup> S GPCR assay	GPCR-mediated guanine nucleotide exchange can be monitored in this assay	500 assays	RPNQ0210

To speak with your local representative, please visit [www.revivity.com/contactus](http://www.revivity.com/contactus).

revvity