

WAVEchips

A sensor chip for every need



EXPERIENCE A NEW
LEVEL OF FLEXIBILITY
IN ASSAY DEVELOPMENT

ROBUST MICROFLUIDICS FOR A BROAD RANGE OF APPLICATIONS



AT CREOPTIX, WE STRIVE FOR FLEXIBILITY AND HIGH PERFORMANCE

The Creoptix WAVEchips offer fast and synchronized sample injections and stability for long off-rates, paired with the broadest sample compatibility and the most user-friendly handling on the market. With no-clog microfluidics and revolutionary sensor technology in one single disposable cartridge, we offer a broad range of sensor chips to best fit your application needs.

FLEXIBILITY

Explore any biomolecular interaction with a broad range of sensor chip surfaces for an all-in-one solution.

COMPATIBILITY

Perform your experiments in biofluids (including serum and plasma), crude samples and uncommon solvents thanks to our no-clog microfluidics design for reduced maintenance costs.

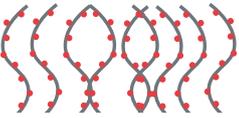
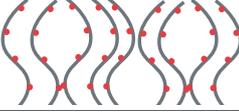
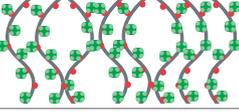
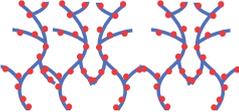
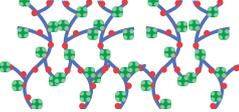
EFFECTIVENESS

Exchange your WAVEchips to have fresh, clean microfluidics within minutes, without the need for special technical support.

PERFORMANCE

Enjoy the fastest and most synchronized sample transitions thanks to our parallel flow channels for high-quality kinetics.

WAVEchips QUICKGUIDE

WAVEchip	Special Characteristics	Matrix		Immobilization Modalities		
		PC	CMD	COV	BIOTIN	HIS
PCH 	Thick hydrogel	✓		✓		
PCP 	Quasi-planar	✓		✓		
PCL 	Thick hydrogel with reduced charges	✓		✓ ₁		
PCH-STA 	Streptavidin-coated	✓			✓	
PCP-STA 	Quasi-planar, streptavidin-coated	✓			✓	
PCH-NTA 	NTA-functionalized	✓				✓
PCP-NTA 	Quasi-planar, NTA-functionalized	✓				✓
PCP-PAG 	Protein A/G-functionalized	✓				
PCP-LIP 	Quasi-planar, lipid anchors	✓				
PCZ 	Zwitterionic polymer	✓		✓		
DXH Upon request 	Thick hydrogel		✓	✓		
DXH-STA Upon request 	Streptavidin-coated		✓		✓	

Legend: polycarboxylate (PC), carboxymethyl dextran (CMD), covalent functionalization through -NH₂, -SH, -CHO, -OH and -COOH (COV), streptavidin/biotin coupling (biotin), His-tag/NTA-coupling (His)



Capacity	Capacity		Suggested Applications	Great for*
	OTHER	HIGH		
		▲	Large ligand-to-analyte molecular weight ratio. General purpose.	   
		▼	Large ligands and/or analytes such as proteo-/liposomes, viruses, VLPs.	  
		▼	Complex matrices such as serum, culture supernatant.	
		▲	Biotinylated ligands. General purpose.	   
		▼	Large biotinylated ligands and/or analytes such as proteo-/liposomes, viruses, VLPs.	   
		▲	His-tagged ligands. General purpose.	  
		▼	Large His-tagged-ligands and/or analytes such as proteo-/liposomes, viruses, VLPs.	 
IgG		▼	Antibody (IgG) ligands.	
Lipid		▼	Hydrophobic ligands such as liposomes, membrane vesicles or fragments. Compatible with large ligands and/or analytes.	 
		▼	Acidic protein and/or negatively charged ligands. Recommended for positively charged analytes. High non-fouling properties.	
		▲	General purpose.	   
		▲	Biotinylated ligands. General purpose.	   

*Applications:

 Small molecules/FBDS

 Serology

 Membrane proteins

 Biologics

Place your order through orders@creoptix.com

For specific questions on our applications, contact us through support@creoptix.com

4PCP
1841A006

creoptix
sensors

EXPLORE NO LIMITS DISCOVER NEW APPLICATIONS

SMALL OR LARGE

From fragment screening to protein, antibody and membrane protein characterization, we have a sensor to suit your needs. Explore a wide range of immobilization modalities and conditions.

FAST OR SLOW

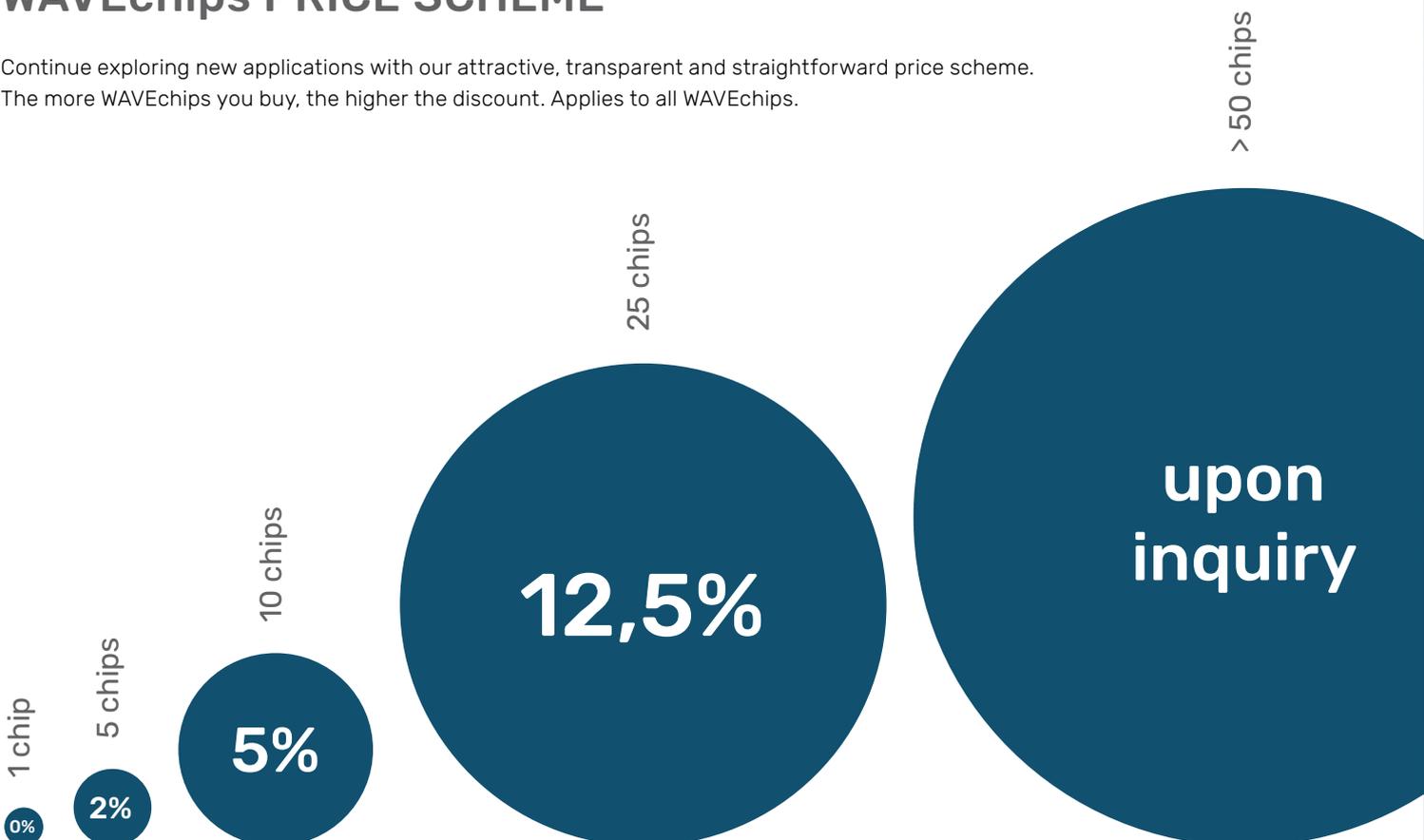
From tight to weak binders, fast to slow off-rates, our microfluidics and surfaces combine long-term stability with highly synchronized and extremely fast transitions from buffer-to-sample and sample-to-buffer.

SIMPLE OR COMPLEX

For extraordinary assay flexibility, our sensors are fully compatible with the toughest samples, including 100% serum or plasma, cell lysates, crude membrane protein preparations, large drug targets, virus-like particles (VLPs), liposomes, and more.

WAVEchips PRICE SCHEME

Continue exploring new applications with our attractive, transparent and straightforward price scheme. The more WAVEchips you buy, the higher the discount. Applies to all WAVEchips.



ORDERING INFORMATION

Choose the most suitable sensor chip for your application
All WAVEchips are suitable for both the WAVE and WAVEdelta systems

POLYCARBOXYLATE-LAYERED SURFACES

WAVECHIP 4PCH SERIES: sensors functionalized with a high capacity polycarboxylate layer

ORDER NO.	Description
4PCH	Sensor chip suited for immobilization through amine coupling
4PCH-STA	Sensor chip with pre-immobilized streptavidin, ideal for capturing biotinylated molecules
4PCH-NTA	Sensor chip with pre-immobilized NTA, ideal for capturing His-tagged molecules

WAVECHIP 4PCP SERIES: sensors functionalized with a thin polycarboxylate layer

ORDER NO.	Description
4PCP	Sensor chip suited for immobilization through amine coupling
4PCP-STA	Sensor chip with pre-immobilized streptavidin, ideal for capturing biotinylated molecules
4PCP-NTA	Sensor chip with pre-immobilized NTA, ideal for capturing His-tagged molecules
4PCP-LIP*	Sensor chip with pre-immobilized lipophilic groups for capturing lipid-based particles or hydrophobic ligands
4PCP-PAG	Sensor chip with pre-immobilized proteinA/G fusion for capturing the broadest range of IgG and FC-tagged proteins

WAVECHIP ALTERNATIVE SERIES: sensors functionalized with special polycarboxylate layers

ORDER NO.	Description
4PCZ	Low fouling sensor functionalized with a zwitterionic polymer bearing carboxylic acids and tertiary amines in similar densities. This sensor chip is suited for immobilization through amine coupling. Ideal for coupling proteins with low isoelectric points or polyanionic ligands
4PCL	Low capacity sensor functionalized with a thick polycarboxylate layer with a lower amount of negative charges (approx. 25%). This sensor chip is suited for immobilization through amine coupling. Requires sulfo-NHS for activation

CARBOXYMETHYL-DEXTRAN-LAYERED SURFACES

WAVECHIP 4DXH SERIES: sensors functionalized with a carboxymethyl-dextran layer

ORDER NO.	Description
4DXH*	High capacity, sensor chip suited for immobilization through amine coupling
4DXH-STA*	High capacity, sensor chip with pre-immobilized streptavidin, ideal for capturing biotinylated molecules

*Available upon request. Longer delivery times may apply
Place your order through orders@creoptix.com

ROBUST MICROFLUIDICS IN A SINGLE DISPOSABLE CARTRIDGE

To stay up-to-date with our latest developments,
visit www.creoptix.com

Place your order through orders@creoptix.com

WAVE⁵



LABEL-FREE DATA
LIKE YOU'VE NEVER SEEN BEFORE

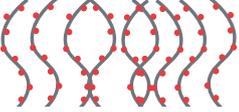
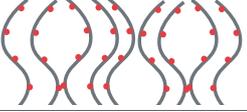
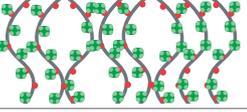
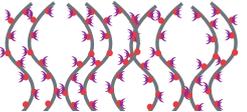
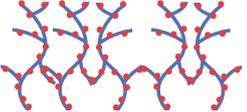
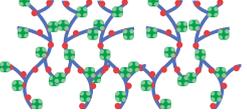


Creoptix AG | Zugerstrasse 76 | 8820 Wädenswil | Switzerland (HQ)
Creoptix Inc | c/o The Brighton Company | 111 South Bedford Street, Ste. 108 | MA-01803 Burlington | USA
www.creoptix.com | +41 44 533 26 60

©2020 Creoptix Sensors / Creoptix. WAVEchips, WAVEcore and waveRAPID are trademarks of CreoptixAG. All goods and support services sold by Creoptix are subject to our terms and conditions. Please visit our website or contact your local Creoptix representative for the most up-to-date information.

KEEPING KINETICS REAL

WAVEchips QUICKGUIDE

WAVEchip	Special Characteristics	Matrix	
		PC	CMD
PCH 	Thick hydrogel	✓	
PCP 	Quasi-planar	✓	
PCL 	Thick hydrogel with reduced charges	✓	
PCH-STA 	Streptavidin-coated	✓	
PCP-STA 	Quasi-planar, streptavidin-coated	✓	
PCH-NTA 	NTA-functionalized	✓	
PCP-NTA 	Quasi-planar, NTA-functionalized	✓	
PCP-PAG 	Protein A/G-functionalized	✓	
PCP-LIP 	Quasi-planar, lipid anchors	✓	
PCZ 	Zwitterionic polymer	✓	
DXH Upon request 	Thick hydrogel		✓
DXH-STA Upon request 	Streptavidin-coated		✓

Immobilization Modality				Capacity		Suggested Applications	Great for*
COV	BIOTIN	HIS	OTHER	HIGH	LOW		
✓				▲		Large ligand-to-analyte molecular weight ratio. General purpose.	
✓					▼	Large ligands and/or analytes such as proteo-/liposomes, viruses, VLPs.	
✓ ¹					▼	Complex matrices such as serum, culture supernatant.	
	✓			▲		Biotinylated ligands. General purpose.	
	✓				▼	Large biotinylated ligands and/or analytes such as proteo-/liposomes, viruses, VLPs.	
		✓		▲		His-tagged ligands. General purpose.	
		✓			▼	Large His-tagged-ligands and/or analytes such as proteo-/liposomes, viruses, VLPs.	
			IgG		▼	Antibody (IgG) ligands.	
			Lipid		▼	Hydrophobic ligands such as liposomes, membrane vesicles or fragments. Compatible with large ligands and/or analytes.	
					▼	Acidic protein and/or negatively charged ligands. Recommended for positively charged analytes. High non-fouling properties.	
✓				▲		General purpose.	
	✓			▲		Biotinylated ligands. General purpose.	



ORDERING INFORMATION

WAVEchips suitable for all WAVE systems

WAVEchip	Quantity	Order No.	Price (USD)
WAVEchip 4PCH	Pack of 1	4PCH-01	295
	Pack of 3	4PCH-03	825
	Pack of 10	4PCH-10	2'595
WAVEchip 4PCP	Pack of 1	4PCP-01	295
	Pack of 3	4PCP-03	825
	Pack of 10	4PCP-10	2'595
WAVEchip 4PCL	Pack of 1	4PCL-01	295
	Pack of 3	4PCL-03	825
	Pack of 10	4PCL-10	2'595
WAVEchip 4PCH-STA	Pack of 1	4PCH-STA-01	395
	Pack of 3	4PCH-STA-03	1'125
	Pack of 10	4PCH-STA-10	3'525
WAVEchip 4PCP-STA	Pack of 1	4PCP-STA-01	395
	Pack of 3	4PCP-STA-03	1'125
	Pack of 10	4PCP-STA-10	3'525
WAVEchip 4PCH-NTA	Pack of 1	4PCH-NTA-01	395
	Pack of 3	4PCH-NTA-03	1'125
	Pack of 10	4PCH-NTA-10	3'525
WAVEchip 4PCP-NTA	Pack of 1	4PCP-NTA-01	395
	Pack of 3	4PCP-NTA-03	1'125
	Pack of 10	4PCP-NTA-10	3'525
WAVEchip 4PCP-PAG	Pack of 1	4PCP-PAG-01	395
	Pack of 3	4PCP-PAG-03	1'125
	Pack of 10	4PCP-PAG-10	3'525
WAVEchip 4PCP-LIP	Pack of 1	4PCP-LIP-01	395
	Pack of 3	4PCP-LIP-03	1'125
	Pack of 10	4PCP-LIP-10	3'525
WAVEchip 4PCZ	Pack of 1	4PCZ-01	395
	Pack of 3	4PCZ-03	1'125
	Pack of 10	4PCZ-10	3'525
WAVEchip 4DXH	Pack of 1	4DXH-01	295
	Pack of 3	4DXH-03	825
	Pack of 10	4DXH-10	2'595
WAVEchip 4DXH-STA	Pack of 1	4DXH-STA-01	395
	Pack of 3	4DXH-STA-03	1'125
	Pack of 10	4DXH-STA-10	3'525



ORDERING INFORMATION

WAVEchips suitable for all WAVE systems

WAVEchip	Quantity	Order No.	Price (EUR)
WAVEchip 4PCH	Pack of 1	4PCH-01	295
	Pack of 3	4PCH-03	825
	Pack of 10	4PCH-10	2'595
WAVEchip 4PCP	Pack of 1	4PCP-01	295
	Pack of 3	4PCP-03	825
	Pack of 10	4PCP-10	2'595
WAVEchip 4PCL	Pack of 1	4PCL-01	295
	Pack of 3	4PCL-03	825
	Pack of 10	4PCL-10	2'595
WAVEchip 4PCH-STA	Pack of 1	4PCH-STA-01	395
	Pack of 3	4PCH-STA-03	1'125
	Pack of 10	4PCH-STA-10	3'525
WAVEchip 4PCP-STA	Pack of 1	4PCP-STA-01	395
	Pack of 3	4PCP-STA-03	1'125
	Pack of 10	4PCP-STA-10	3'525
WAVEchip 4PCH-NTA	Pack of 1	4PCH-NTA-01	395
	Pack of 3	4PCH-NTA-03	1'125
	Pack of 10	4PCH-NTA-10	3'525
WAVEchip 4PCP-NTA	Pack of 1	4PCP-NTA-01	395
	Pack of 3	4PCP-NTA-03	1'125
	Pack of 10	4PCP-NTA-10	3'525
WAVEchip 4PCP-PAG	Pack of 1	4PCP-PAG-01	395
	Pack of 3	4PCP-PAG-03	1'125
	Pack of 10	4PCP-PAG-10	3'525
WAVEchip 4PCP-LIP	Pack of 1	4PCP-LIP-01	395
	Pack of 3	4PCP-LIP-03	1'125
	Pack of 10	4PCP-LIP-10	3'525
WAVEchip 4PCZ	Pack of 1	4PCZ-01	395
	Pack of 3	4PCZ-03	1'125
	Pack of 10	4PCZ-10	3'525
WAVEchip 4DXH	Pack of 1	4DXH-01	295
	Pack of 3	4DXH-03	825
	Pack of 10	4DXH-10	2'595
WAVEchip 4DXH-STA	Pack of 1	4DXH-STA-01	395
	Pack of 3	4DXH-STA-03	1'125
	Pack of 10	4DXH-STA-10	3'525



ORDERING INFORMATION

WAVEchips suitable for all WAVE systems

WAVEchip	Quantity	Order No.
WAVEchip 4PCH	Pack of 1 Pack of 3 Pack of 10	4PCH-01 4PCH-03 4PCH-10
WAVEchip 4PCP	Pack of 1 Pack of 3 Pack of 10	4PCP-01 4PCP-03 4PCP-10
WAVEchip 4PCL	Pack of 1 Pack of 3 Pack of 10	4PCL-01 4PCL-03 4PCL-10
WAVEchip 4PCH-STA	Pack of 1 Pack of 3 Pack of 10	4PCH-STA-01 4PCH-STA-03 4PCH-STA-10
WAVEchip 4PCP-STA	Pack of 1 Pack of 3 Pack of 10	4PCP-STA-01 4PCP-STA-03 4PCP-STA-10
WAVEchip 4PCH-NTA	Pack of 1 Pack of 3 Pack of 10	4PCH-NTA-01 4PCH-NTA-03 4PCH-NTA-10
WAVEchip 4PCP-NTA	Pack of 1 Pack of 3 Pack of 10	4PCP-NTA-01 4PCP-NTA-03 4PCP-NTA-10
WAVEchip 4PCP-PAG	Pack of 1 Pack of 3 Pack of 10	4PCP-PAG-01 4PCP-PAG-03 4PCP-PAG-10
WAVEchip 4PCP-LIP	Pack of 1 Pack of 3 Pack of 10	4PCP-LIP-01 4PCP-LIP-03 4PCP-LIP-10
WAVEchip 4PCZ	Pack of 1 Pack of 3 Pack of 10	4PCZ-01 4PCZ-03 4PCZ-10
WAVEchip 4DXH	Pack of 1 Pack of 3 Pack of 10	4DXH-01 4DXH-03 4DXH-10
WAVEchip 4DXH-STA	Pack of 1 Pack of 3 Pack of 10	4DXH-STA-01 4DXH-STA-03 4DXH-STA-10



Creoptix AG | Zugerstrasse 76 | 8820 Wädenswil | Switzerland (HQ)

Creoptix Inc | c/o The Brighton Company | 111 South Bedford Street, Ste. 108 | MA-01803 Burlington | USA

www.creoptix.com | +41 44 533 26 60