

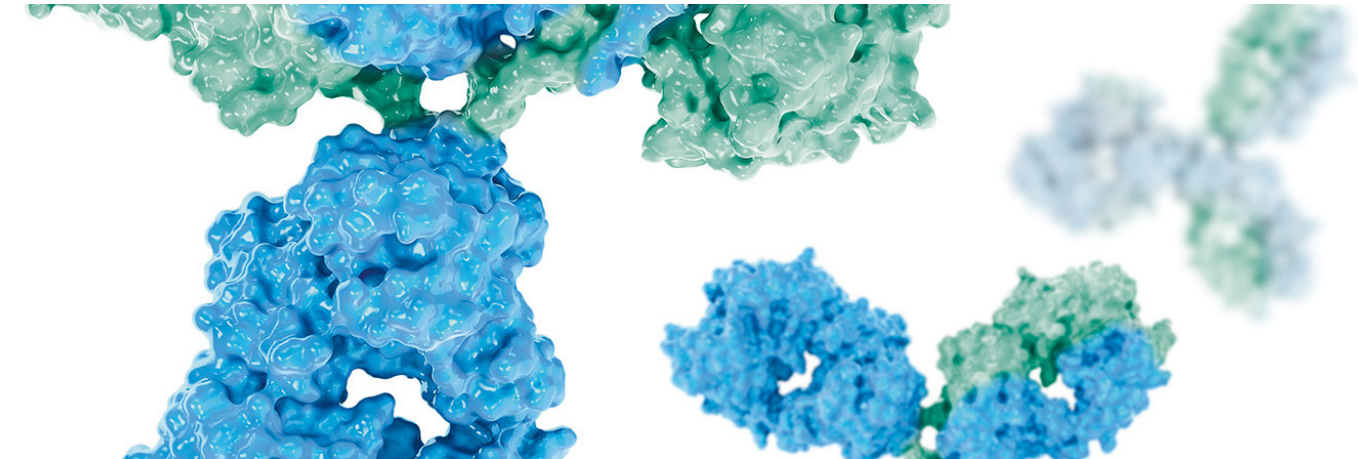
The background of the entire page is a 3D molecular model of a protein complex, likely a tetramer, rendered in shades of blue and yellow. The protein is shown in a surface representation, highlighting its complex, multi-subunit structure. The background is a dark blue gradient with some light blue bokeh effects.

Invitation  
MicroCal Experience Day  
15 September 2022

# Program

09.30 - 10.00	Registration and coffee
10.00 - 10.15	Welcome - <i>Tom Daneels - Jan Volkaert, Sysmex</i>
10.15 - 10.45	MicroCal PEAQ-ITC: the gold standard for measuring binding affinities - <i>Kris Rogiers, Sysmex</i>
10.45 - 11.15	Analysis of molecular glues with ITC <i>Loes Stevers, Ambagon Therapeutics</i>
<b>BREAK</b>	
11.30 - 12.00	Structural and thermodynamic insights into protein-DNA interactions using ITC in combination with other biophysical techniques - <i>Maruša Prolic Kalinsek, VUB</i>
12.00 - 12.15	ITC hands-on session - <i>Kris Rogiers, Sysmex</i>
<b>LUNCH</b>	
13.00 - 13.30	MicroCal PEAQ-DSC: the gold standard for protein stability characterization <i>Peter Gimeson, Malvern Panalytical</i>
13.30 - 14.00	Delving deeper into AAV attributes using DSC - <i>Peter Gimeson, Malvern Panalytical</i>
<b>BREAK</b>	
14.15 - 14.30	DSC hands-on session - <i>Peter Gimeson, Malvern Panalytical</i>
14.30 - 15.00	Pushing the Boundaries in Biomolecular Interaction Analysis with the Creoptix WAVEsystem. From Biologics to Fragment-Based Drug Discover - <i>Application Scientist, Creoptix/Malvern Panalytical</i>
15.00 - 15.30	Protein Interaction for precision medicines development using Creoptix WAVEsystem <i>Ellen Mattaar, Byondis</i>
<b>BREAK</b>	
15.45 - 16.15	Binding kinetics with Creoptix WAVEsystem for innovative drug development <i>David Moreno Delgado, Galapagos</i>
16.15 - 16.30	WAVE system hands-on session - <i>Application Scientist, Creoptix/Malvern Panalytical</i>
16.30 - 16.45	Q&A
16.45	Closing

# MICROCAL EXPERIENCE DAY for all ins and outs on microcalorimetry



We would like to invite you to our MicroCal Experience Day on the 15th of September at our office in Etten-Leur (The Netherlands).

Both MicroCal techniques, ITC (Isothermal Titration Calorimetry) and DSC (Differential Scanning Calorimetry), will be discussed with tips and tricks through presentations, guest lectures, and hands-on sessions.

In addition, the new GCI (Grating-Coupled Interferometry) technology from Creoptix will also be presented.

We look forward to seeing you on September 15th!

Kind regards,  
Your Sysmex I&R Team

Register here:





---

**Sysmex Nederland B.V.**

Ecustraet 11, 4879 NP Etten-Leur, The Netherlands · Phone +31 76 5086000 · Fax +31 76 5086086 · info@sysmex.nl · [www.sysmex.nl](http://www.sysmex.nl)

**Sysmex Belgium N.V.**

Park Rozendal, Building A, Terhulpsessesteenweg 6a, 1560 Hoeilaart, Belgium · Phone +32 2 7697474 · Fax +32 2 7697499 · info@sysmex.be · [www.sysmex.be](http://www.sysmex.be)