

Biologics Drug Development

Reliable solutions to support you at every step.



Discovery and Early Development

● Hit-to-lead: Hit validation

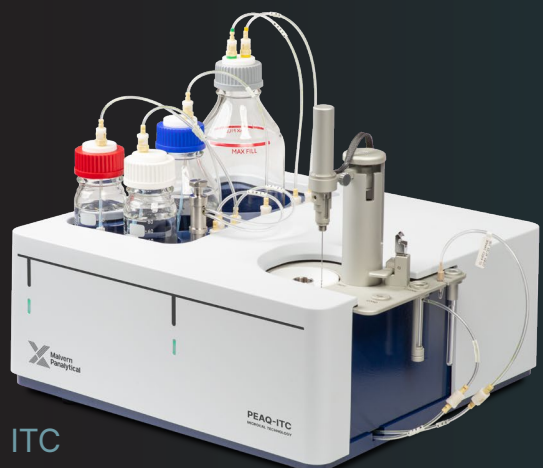
 **Binding affinity & kinetics:**
GCI, ITC

 **Particle size:**
DLS

 **Thermal stability & higher order structure:**
DSC



GCI
WAVEsystem



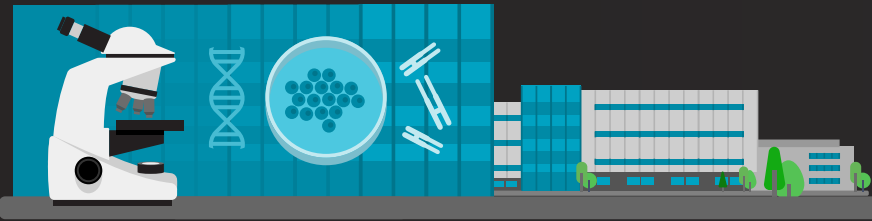
ITC
MicroCal PEAQ ITC



DLS
Zetasizer Advance



DSC
MicroCal PEAQ-DSC



Preclinical Development

● Lead/candidate optimization

 **Particle size, shape & concentration:**
DLS, NTA

 **Thermal stability & higher order structure:**
DSC



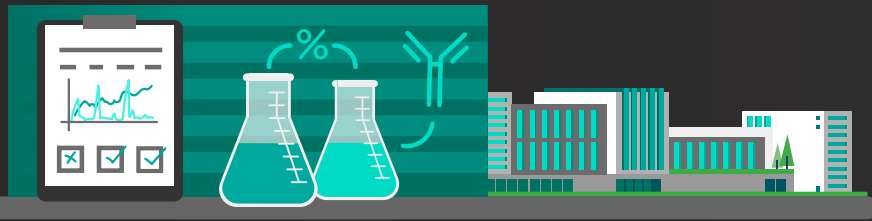
NTA
NanoSight Pro



DLS
Zetasizer Advance



DSC
MicroCal PEAQ-DSC



Clinical & Scale-up

● Formulation & process:
Development & optimization

 **Colloidal stability:**
ELS

 **Thermal stability & higher order structure:**
DSC

 **Particle size, shape & identification (ID):**
DLS, NTA, MDRS



ELS & DLS
Zetasizer Advance



DSC
MicroCal PEAQ-DSC



NTA
NanoSight Pro



MDRS
Morphologi 4-ID



Manufacturing

● QA/QC

 **Thermal stability & higher order structure:**
DSC

 **Particle size & identification (ID):**
DLS, MDRS



DSC
MicroCal PEAQ-DSC



DLS
Zetasizer Advance



MDRS
Morphologi 4-ID