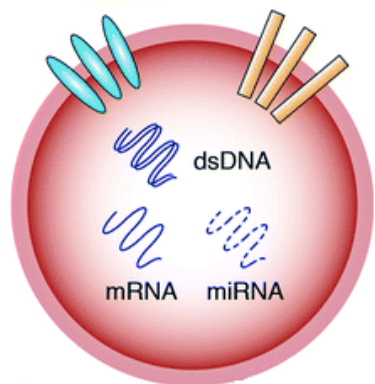


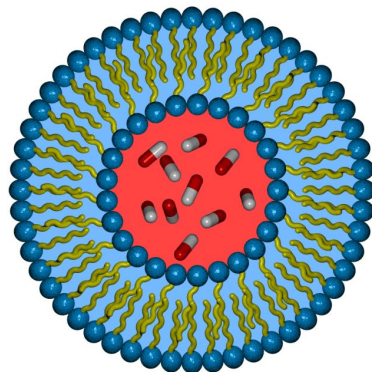
Improved Methods For Fluorescent Labelling And Detection with Nanoparticle Tracking Analysis (NTA)

Aymeric AUDFRAY, PhD
Field Application Scientist, Pharma & Food, France
aymeric.audfray@malvern.com

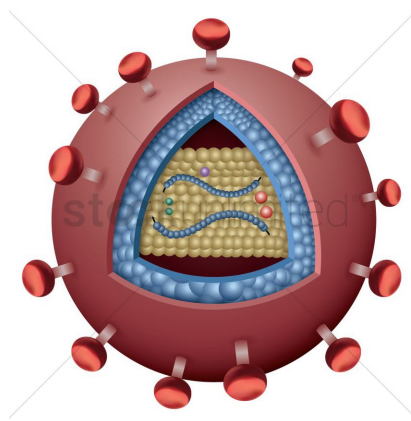
Exosomes



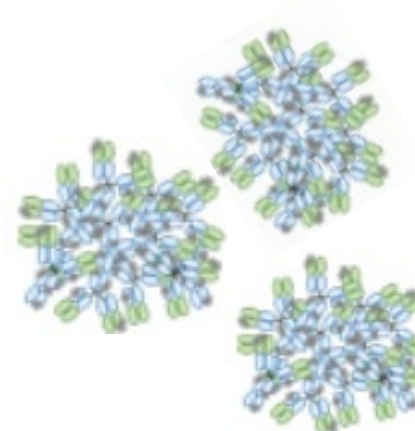
LNPs
(Nanoparticles)



Viruses



Protein Aggregates



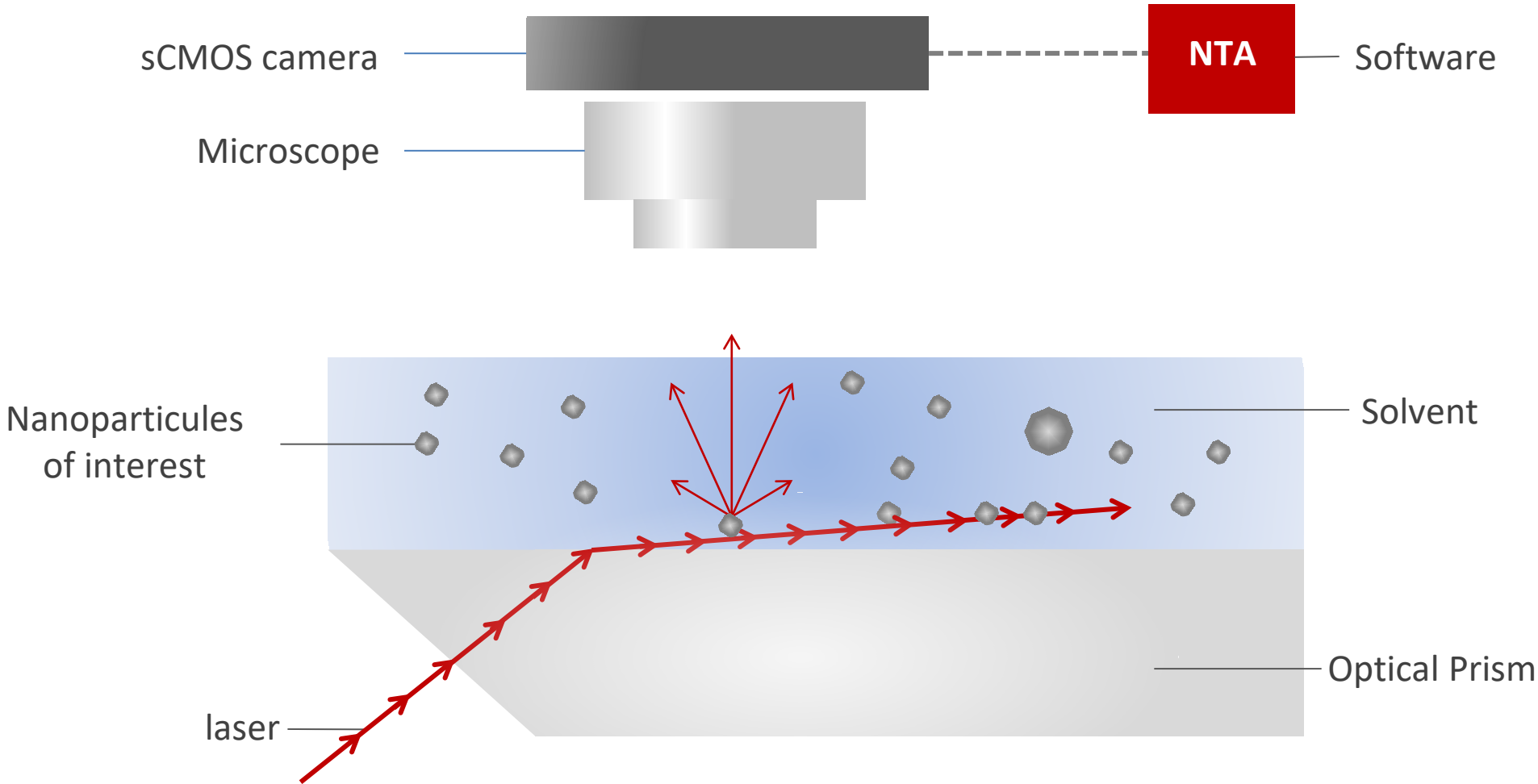
Nanobubbles



Malvern Panalytical



Principle of Nanoparticle Tracking Analysis



Patent:
Bob Carr
(CEO Nanosight 2003)

PCT/GB03/001827

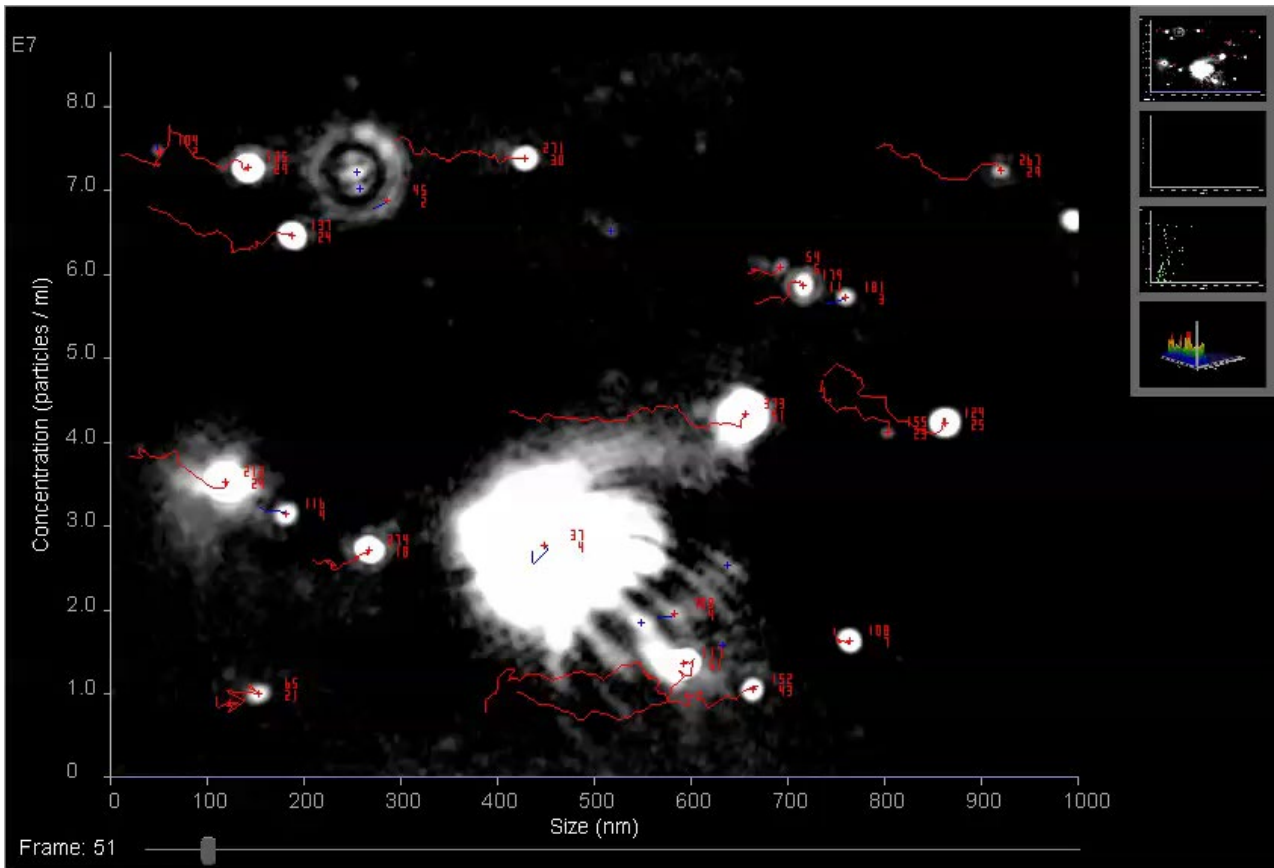


Principle of Nanoparticle Tracking Analysis

Video-captures = raw data

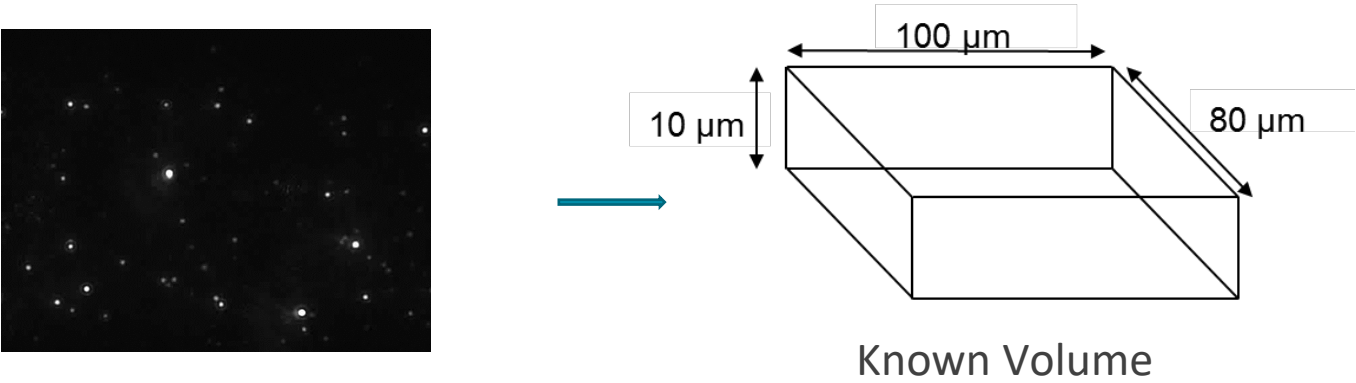


Processing = Tracking of Brownian Motion

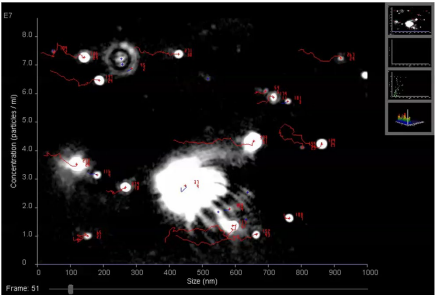


Principle of Nanoparticle Tracking Analysis

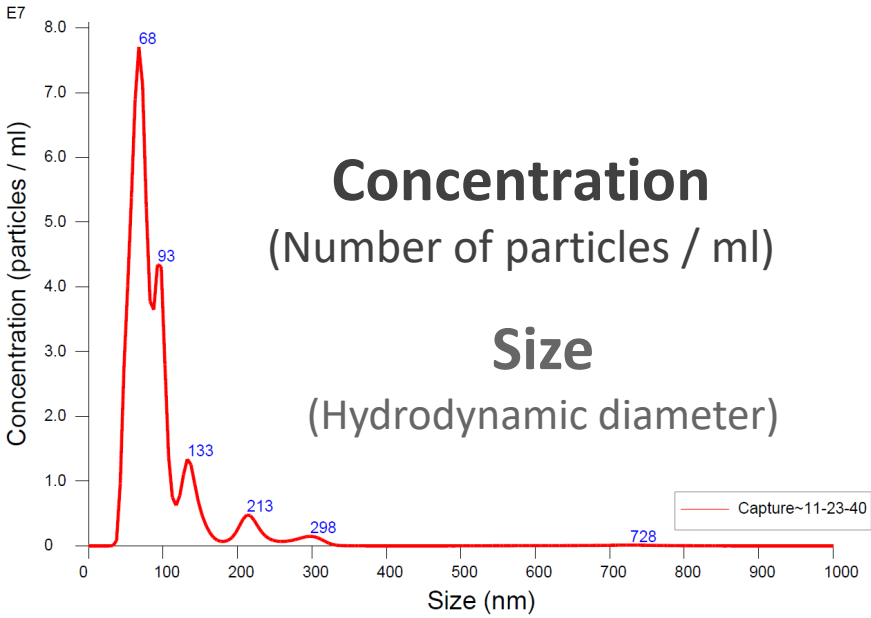
Average number of particles / frame



Tracking of Brownian Motion
= Diffusion Coefficient (m²/s)



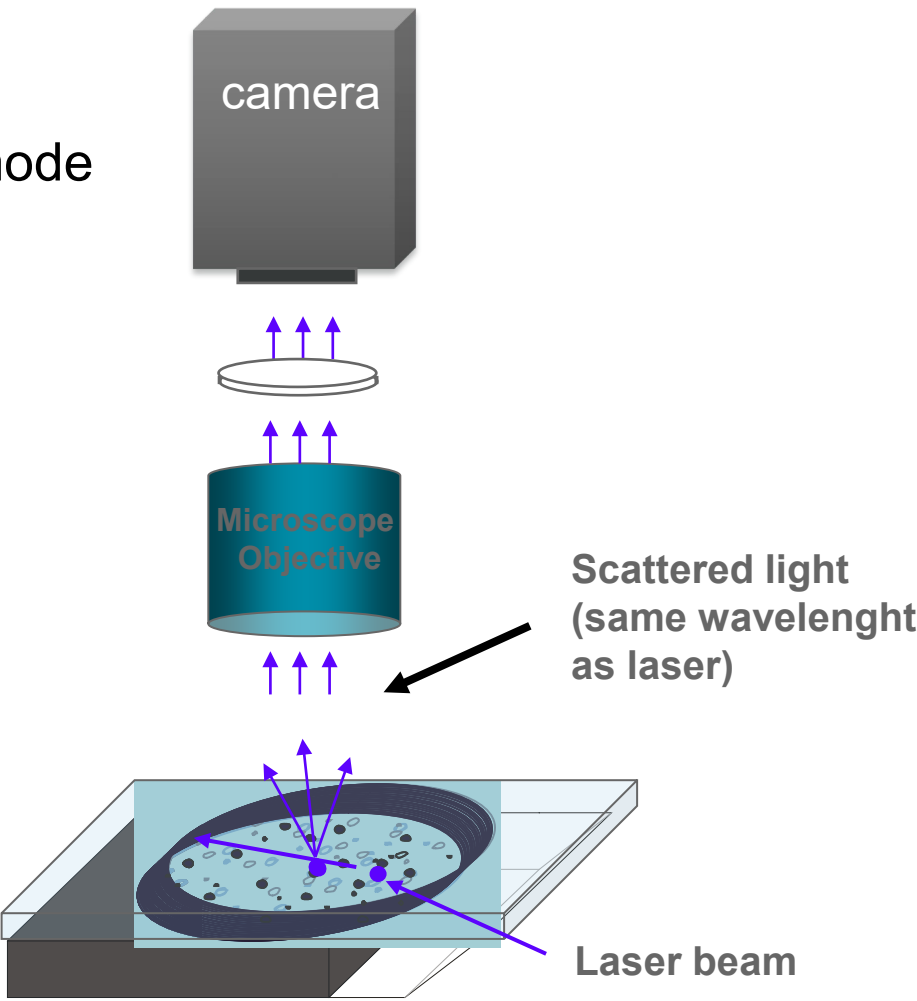
$$D_t = \frac{K_B T}{3\pi\eta d_h}$$



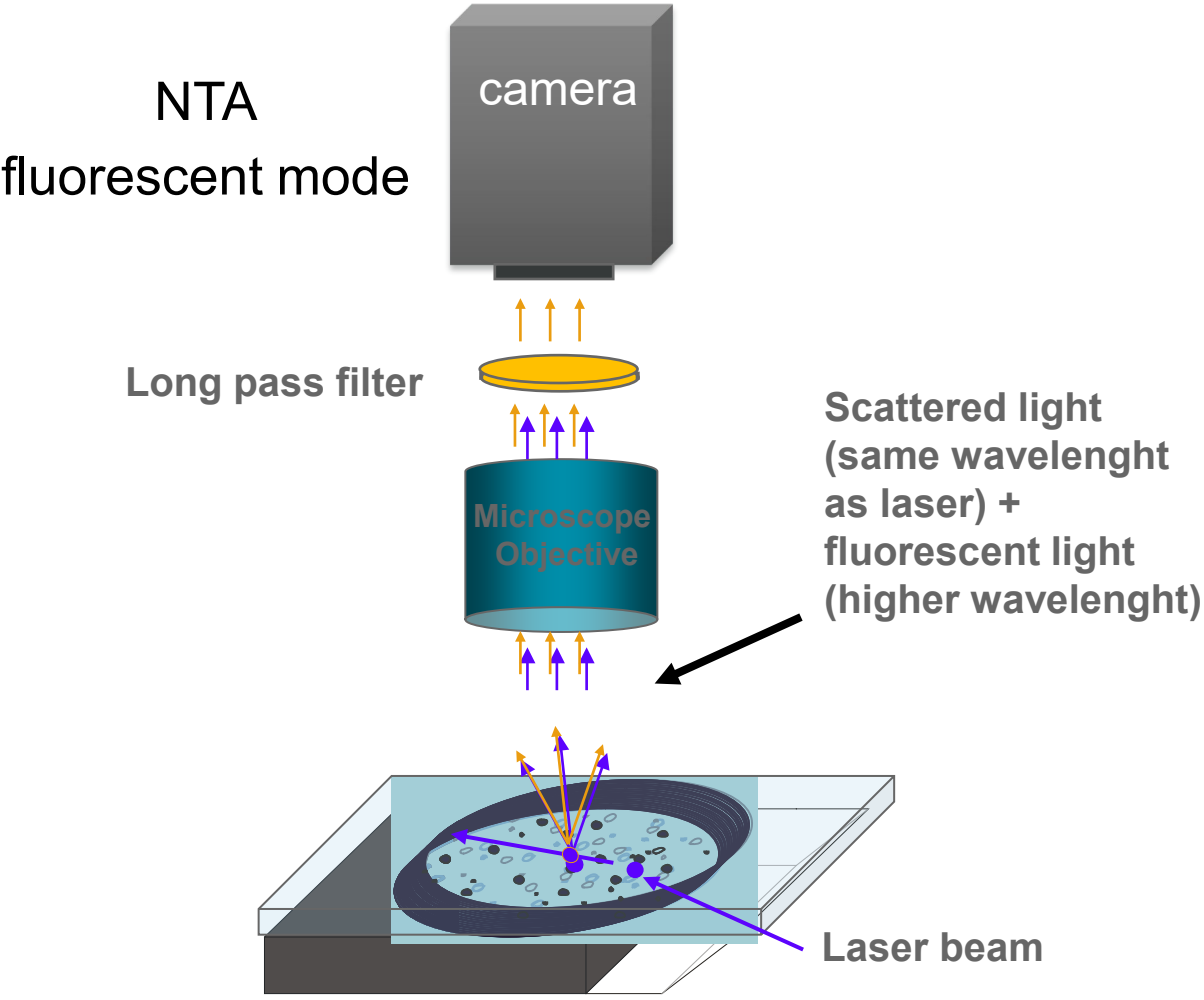
NTA vs Fluo-NTA

Principle

NTA
scatter mode



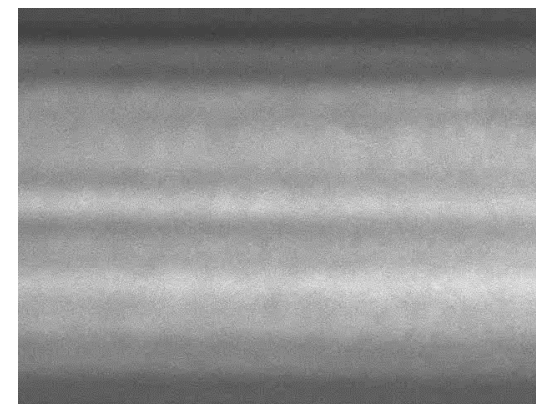
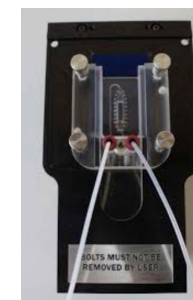
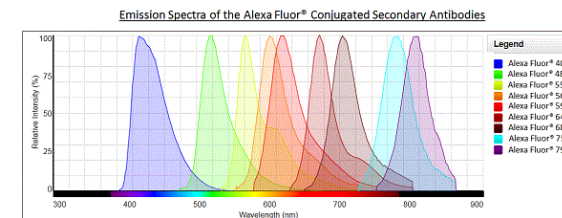
NTA
fluorescent mode



General Tips for Fluorescent Measurements



- Fluo-NTA is always more difficult than scatter mode
 - Samples are more complex (labelling) and settings are more complicated (signal)
- Minimum number of fluorescent dyes by particle:
 - 30-50 for small molecule dyes
 - 3-5 for Quantum Dots
 - Dye excitation maxima must match laser wavelength (405nm, 488nm, 532nm, 642nm)
- Do all operations likely to reduce photobleaching:
 - Trigger cable (present in all instruments now)
 - Use syringe pump
 - Connect inlet tubing to the right port of the LVFC
 - Protect syringe, tubings,... from light.
- Minimize unbound fluorescent dye (generate background noise)
- Check focus setting (different focal plan between scatter and fluorescence)
- Be careful at the dilution (depends on the % of labelling)
- Use advanced camera settings

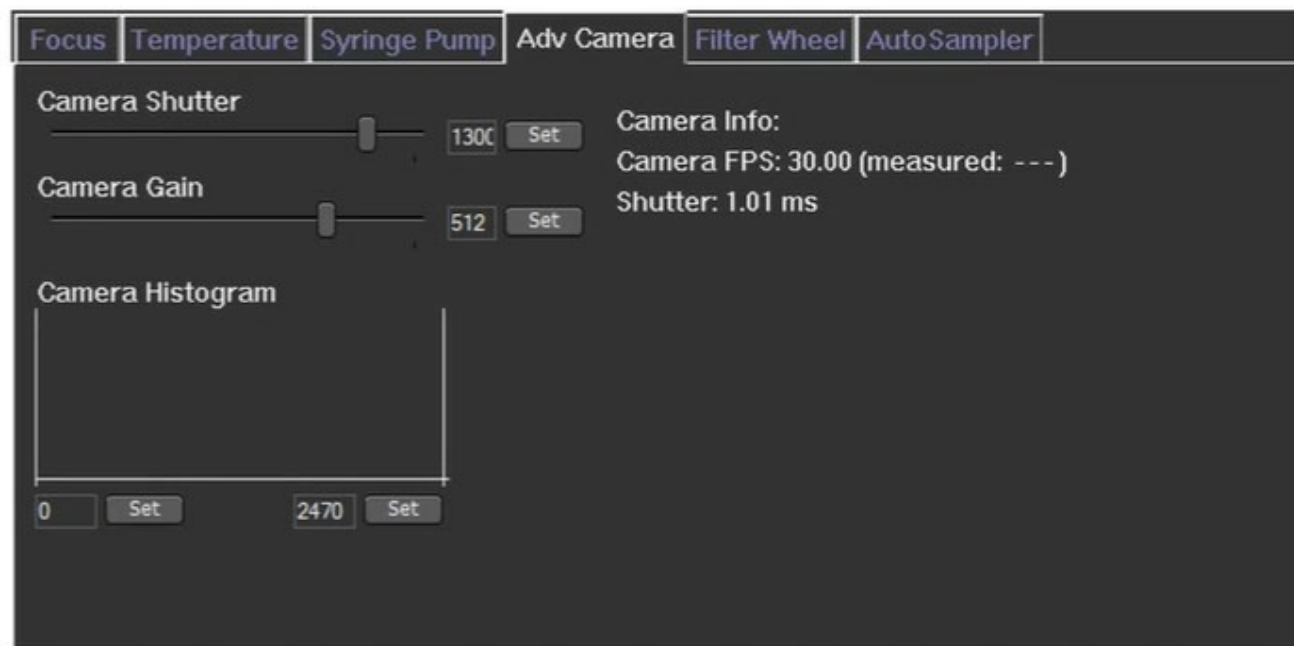


Advanced Camera Settings

Fluo-NTA



- Usually CL is at 16 (maximum) for fluorescent NTA
- Possibility to increase the signal with advanced camera settings
 - Push Camera Shutter/Camera Gain to the right
 - Adjust left (contrast) and right (light) threshold of Camera Histogram



Labelling Strategies

Fluo-NTA



- EVs labelling is the most common application for fluo-NTA
- Membrane Labelling:
 - Exoglow (488nm laser, <https://systembio.com/shop/exoglow-nta-fluorescent-labelling-kit/>)
 - CellMaskOrange (532nm laser)
 - Specific to EVs over protein aggregates, easy labelling with hundreds of dyes by particle
 - Non specific to surface markers
- Antibody labelling
 - Highly specific to surface biomarkers
 - Can be challenging to obtain sufficient signal (30-50 tags per particle)

Fluorescent Labelling of membranes and analysis by Malvern Panalytical Fluo-NTA

Aymeric AUDFRAY, PhD
Field Application Scientist, Pharma & Food, France
aymeric.audfray@malvern.com

NanoSight Characterisation of Fractionated Exosomes (EVs) Labelled with CellMask™ Orange

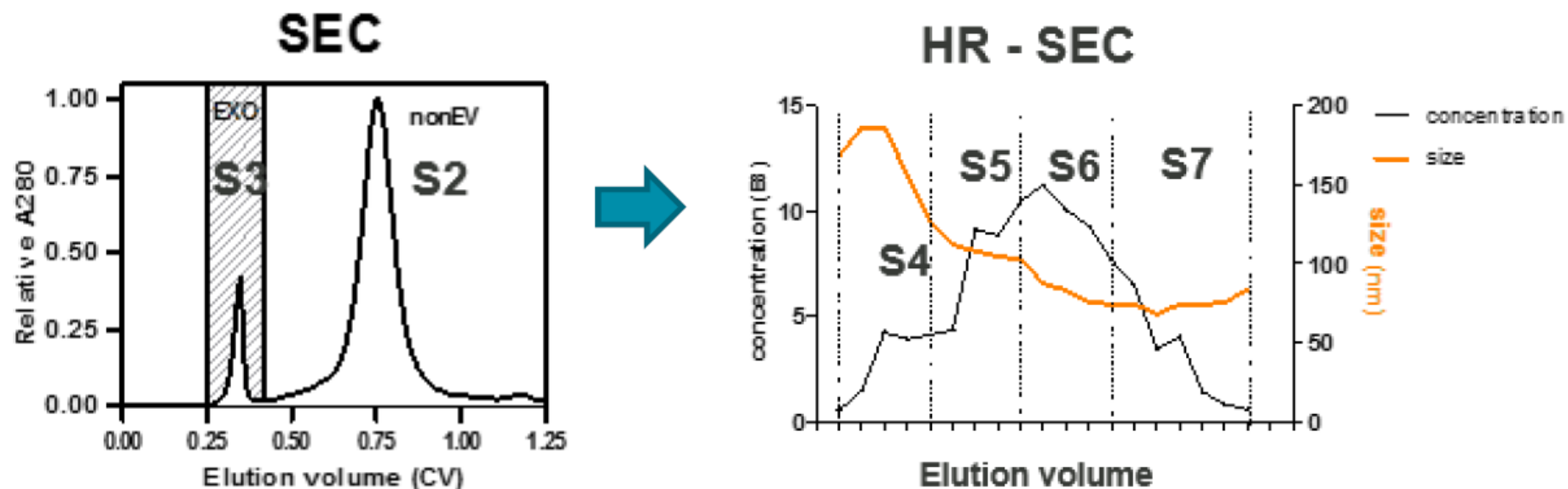
Agnieszka Siupa¹, Eduard Willms², Pauline Carnell¹, Imre Mager²

¹Malvern Panalytical, Grovewood Road, WR14 1XZ, Malvern, Worcestershire, UK

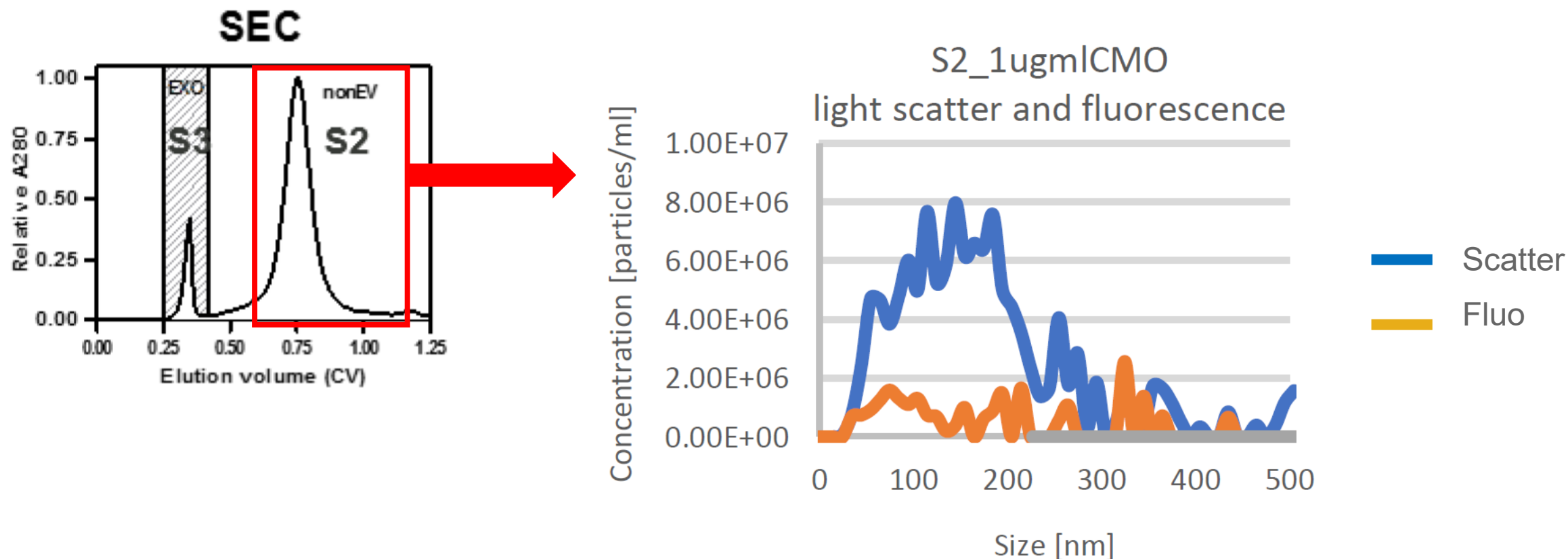
²Department of Physiology, Anatomy and Genetics, Le Gros Clark Building, South Parks Road, Oxford

Purification of EVs

- EVs were isolated from cell culture supernatant of SKOV3 cells
- Differential centrifugation approach
- Size exclusion chromatography (**SEC**) /High resolution size exclusion chromatography (**HR-SEC**)

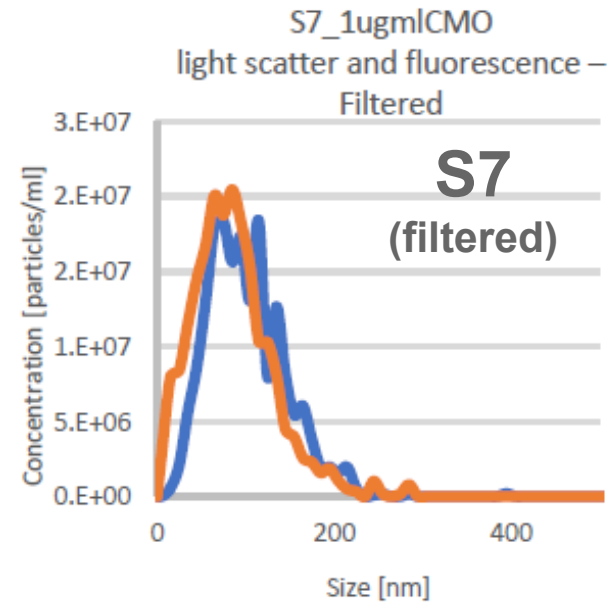
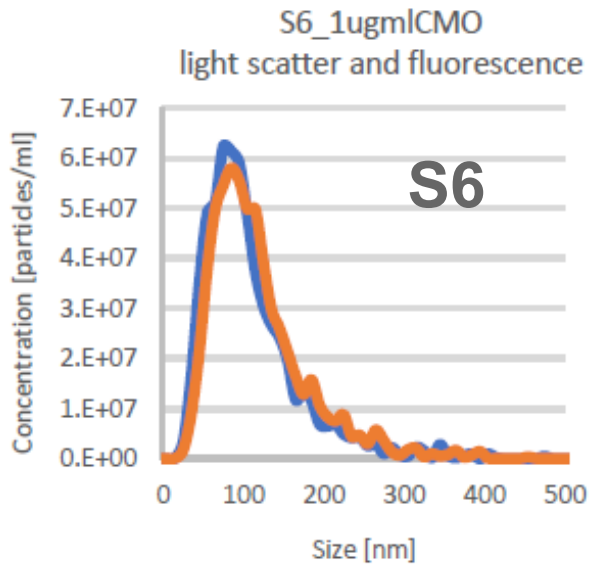
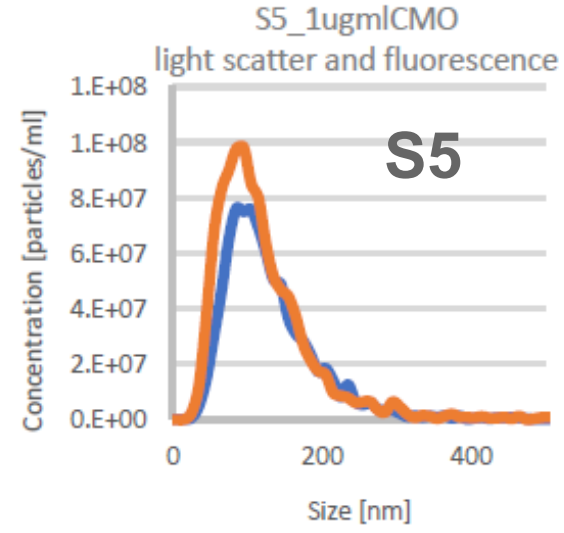
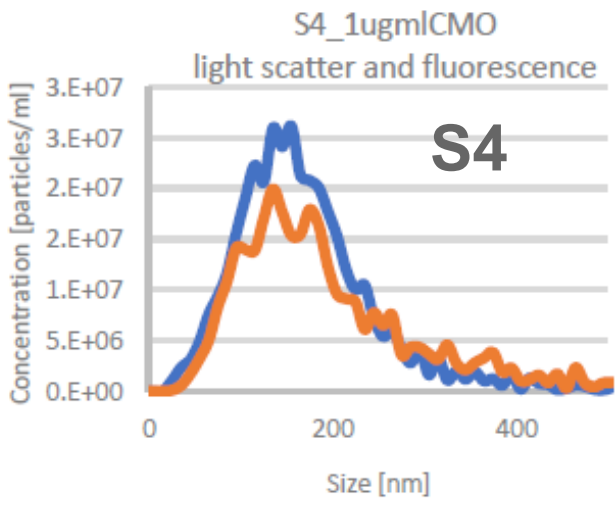
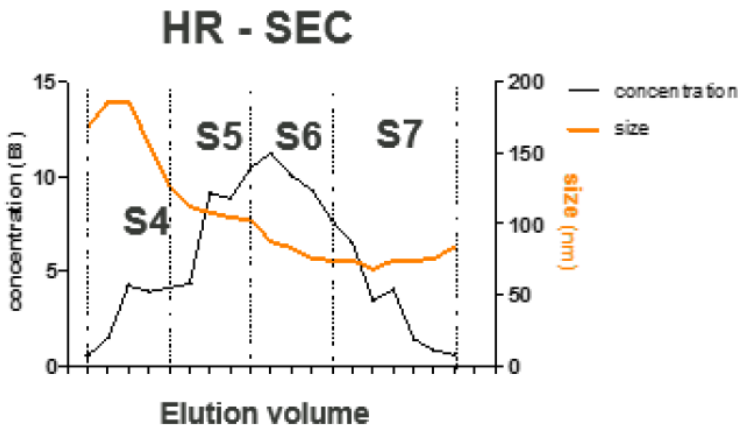


Labelling of EVs with CellMaskOrange® (CMO)



- Fraction S2 contains Nanoparticles that absorbs UV (most likely protein aggregates)
- **No fluorescence signal**
- **No labelling by CMO**

Comparison of scatter and fluo results



— Scatter
— Fluo

Conclusion

- Efficient and simple method to analyse **only** EVs in a samples using Fluo-NTA
- Universal Protocol developped by Malvern Panalytical – **Available on demand**
- Standardization of EVs concentration and size analysis
- Other protocol are being developped with different dyes and wavelenghts


OPEN

Improved methods for fluorescent labeling and detection of single extracellular vesicles using nanoparticle tracking analysis

Received: 24 January 2018

Accepted: 26 July 2019

Published online: 23 August 2019

Kristen E. Thane, Airl M. Davis & Andrew M. Hoffman 

**SCIENTIFIC
REPORTS**
nature research




- EVs extracted from:
 - Canine Mesenchymal Stromal cells (MSCs) of placental origin
 - Human Embryonic Kidney cells ATCC CRL-1573 (HEK-293)
- Isolation using Ultracentrifugation
- Stored at 4°C before immunolabelling (0-4 days)
- NTA Measurement:
 - NS300 with syringe pump
 - Laser 488nm + long-pass filter 500nm
 - SOP: 5x30-60s with flow mode (minimum of 1000 valid tracks)

OPEN Improved methods for fluorescent labeling and detection of single extracellular vesicles using nanoparticle tracking analysis

Received: 24 January 2018

Accepted: 26 July 2019

Published online: 23 August 2019

Kristen E. Thane, Airl M. Davis & Andrew M. Hoffman 

SCIENTIFIC
REPORTS
nature research



- Labelling protocol:
 - Direct:
 - anti-CD9-Qd655
 - anti-CD81-APC
 - Indirect:
 - anti-CD9 + Secondary Ab-Qd655
 - **anti-CD9-biotin + Streptavidin-Qd655**
 - anti-CD81 + Secondary Ab-Qd655
 - anti-CD81-biotin + Streptavidin-Qd655
- 10^{10} EVs + 1 μ g of Anti-CD in 100 μ l of PBS, 4°C Overnight,
- Ultracentrifugation 100 000 g, 2 hours and resuspension in 50-100 μ l
- Anayzed immediatly (after dilution)

OPEN

Improved methods for fluorescent labeling and detection of single extracellular vesicles using nanoparticle tracking analysis

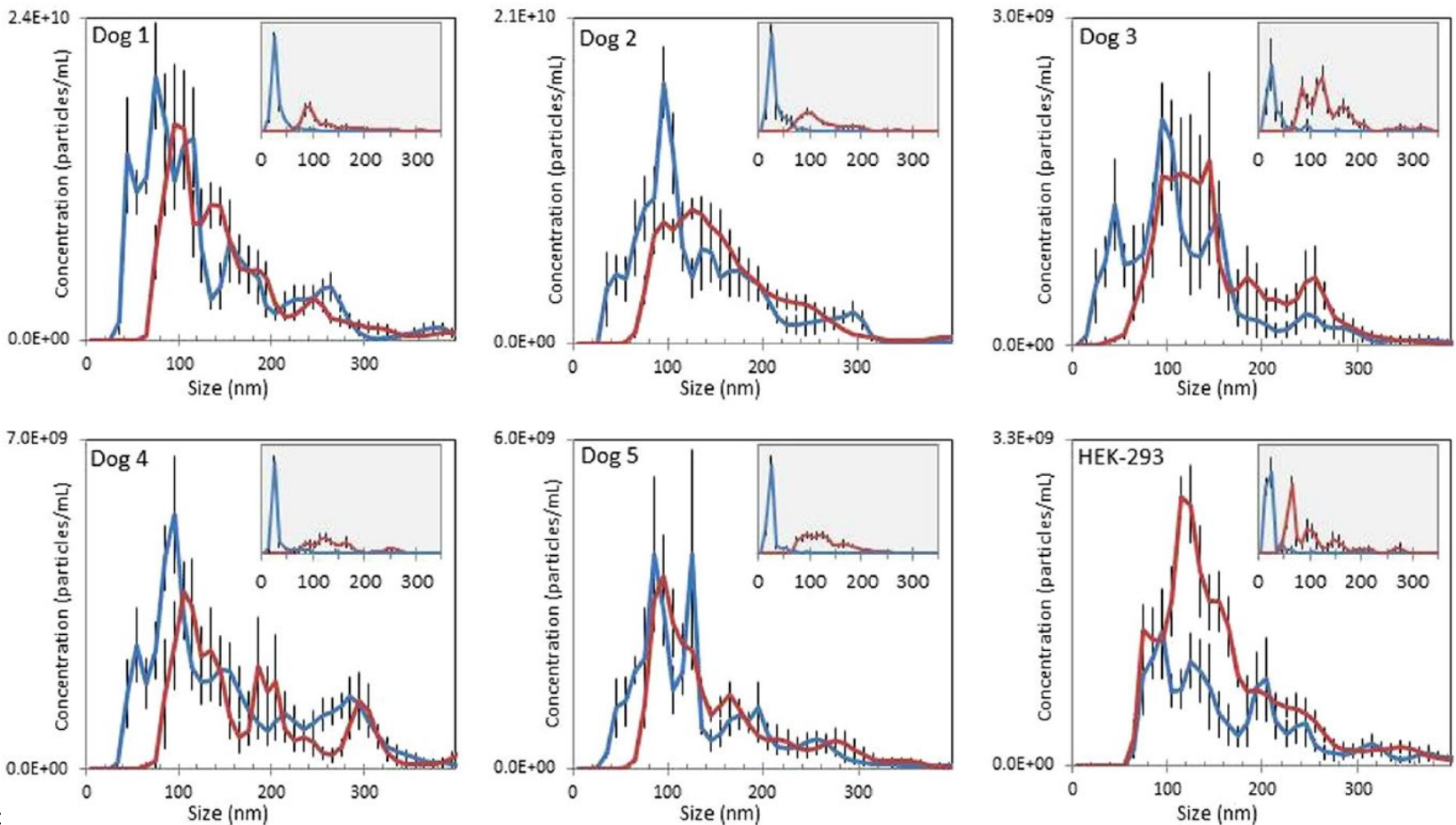
SCIENTIFIC
REPORTS
nature research



Received: 24 January 2018
Accepted: 26 July 2019
Published online: 23 August 2019

Kristen E. Thane, Airlie M. Davis & Andrew M. Hoffman

- anti-CD9-biotin + Streptavidin-Qd655 :



Scatter Mode

Fluo Mode

OPEN

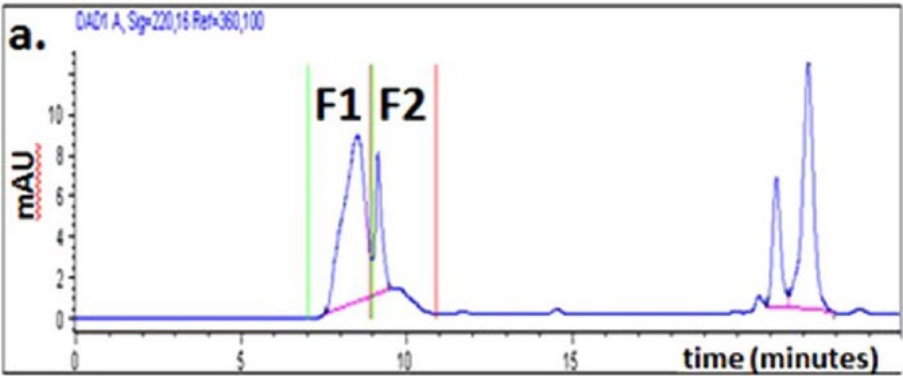
Improved methods for fluorescent labeling and detection of single extracellular vesicles using nanoparticle tracking analysis

Received: 24 January 2018
Accepted: 26 July 2019
Published online: 23 August 2019

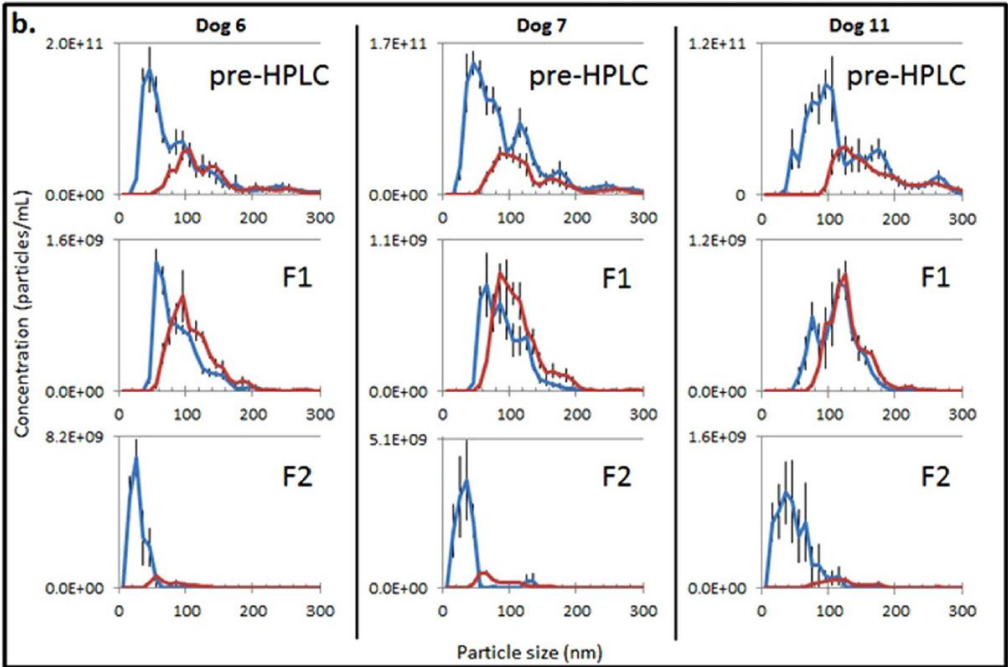
Kristen E. Thane, Airl M. Davis & Andrew M. Hoffman



- Purification of free Streptavidin-Qd655
- 2 methods:
 - Capture of Strep-Qdot using Biotinylated-beads (not shown here)
 - HPLC



Fluo Mode Scatter Mode




OPEN

Improved methods for fluorescent labeling and detection of single extracellular vesicles using nanoparticle tracking analysis

Received: 24 January 2018

Accepted: 26 July 2019

Published online: 23 August 2019

Kristen E. Thane, Airl M. Davis & Andrew M. Hoffman 

**SCIENTIFIC
REPORTS**
nature research



- Conclusion:
 - Development of a method to standardize EVs characterization using Fluo-NTA (protocol and analysis)
 - Use of Quantum dots with 488nm laser
 - Best approach: CD9-biot + Streptavidin-Qd655
 - Develop a method to remove free fluorophore using HPLC or Biotinylated-beads

General Conclusion



- Scatter mode NTA is already used as routine for NPs characterization
- Fluo-NTA can be more challenging (depending on samples)
- Different labelling approaches (membranes, markers,...)
 - CellMaskOrange labelling is universal and available on demand
 - Protocols from bibliography
 - Other specific protocol have to be developped
- Importance of labelling protocol to have sufficient signal/decrease photobleaching
 - One good example: indirect labelling using **Ab-biot + Streptavidin-Qdots**
 - Use of **Qdots** seems very promising
- Support from Malvern Panalytical
- Support from « Concept Life Science » for method development



THANKS FOR YOUR ATTENTION



Labelling of EVs with CellMaskOrange[®] (CMO)



- CellMaskOrange[®] is a dye able to bind specifically to lipid bilayers
- 10µl of 1µg/ml CMO + 10µl EVs (non diluted)
- Incubation 10min at room temperature
- Dilution of samples to optimal concentration range for NTA measurements
- Analyzed on NS300 device
 - 532 nm laser
 - 565 nm fluorescence filter
 - Flow mode
 - Concentration upgrade
- Analysis have been made using both scatter and fluorescence mode