

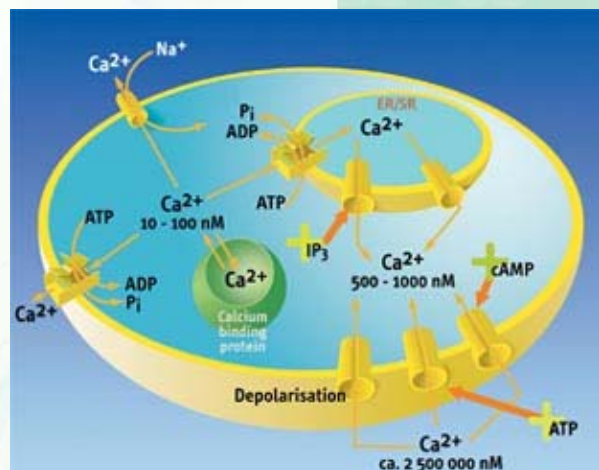
Calcium Monitoring with Mithras LB 940 multimode microplate reader



- up to 4 reagent injectors
- temperature control
- flexible software for operation
- sophisticated kinetics data reduction
- high sensitivity
- ratiometric readout
- fast filter change

Ca²⁺ is an intracellular messenger in many eukaryotic signal transduction pathways. Most Ca²⁺-signalling systems have one thing in common: they generate brief pulses of Ca²⁺, thereby regulating cellular functions.

[Download the brand new Application Note about Calcium Monitoring](#)



[Learn more about Mithras LB 940](#)

[Learn more about Calcium monitoring](#)

With [Interchim](#) - a supplier of high quality research reagents - Berthold Technologies has formed a strong alliance to benefit you as a user.



[Get in touch with Berthold Technologies](#)

www.Berthold.com/Bio