

Microemulsion characterisation using dynamic light scattering



Introduction

Microemulsions have attracted large interest in the pharmaceutical industry as drug delivery systems due to their improved drug solubilisation properties, increased shelf life and ease of preparation. They normally consist of an aqueous phase, an oil phase, a surfactant and a co-surfactant (usually an alcohol). When the concentrations of these components are favorable, they spontaneously emulsify to form a monodisperse, thermodynamically stable, transparent microemulsion.

Many drugs that are insoluble in aqueous media are prepared as pre-concentrates consisting of oil, surfactant and alcohol. This is then diluted with water to form the microemulsion prior to administration to the patient.

Size characterisation of the resulting microemulsion is essential in ensuring safe and efficient dosage. Monitoring of changes in the size distribution can provide valuable information for optimizing the formulation.

Dynamic light scattering (DLS) is a technique for measuring the particle size of colloidal suspensions. In DLS, the sample is illuminated with a laser beam and the intensity of the resulting scattered light produced by the particles fluctuates at a rate that is dependent upon the size of the particles. Analysis of these intensity fluctuations yields the diffusion coefficient of the particles and hence the particle size. Further information on the technique of dynamic light scattering can be found on the Malvern web site.

Experimental

The measurements shown in this application note demonstrate the use of DLS for studying the effect of incorporation of drug into a microemulsion and the time-dependent loss of drug solubility.

Measurements were made on a Zetasizer Nano ZS instrument at 25°C. The instrument contains a 4 mW He-Ne laser operating at a wavelength of 633 nm and incorporates non-invasive backscatter optics (NIBS). The measurements were made at a detection angle of 173° and the measurement position within the cuvette was automatically determined by the software.

The pre-concentrate formulations were diluted in deionised water to form the microemulsion.

Results and Discussion

Figure 1 shows the intensity size distributions obtained from three repeat measurements of the microemulsion. This sample did not contain any drug material. The results show very repeatable intensity size

distributions with z-average diameters of 18.9nm. The z-average diameter is the mean hydrodynamic diameter and is calculated according to the International Standard on dynamic light scattering ISO13321.

Figure 2 shows the intensity size distribution of a microemulsion obtained by dilution of a pre-concentrate formulation in deionised water. This sample contains an oil-soluble drug. The first peak at around 30nm arises from the microemulsion droplets. The peak positioned at 2 microns arises from insoluble drug leaking out of the microemulsion. The presence of the second peak has resulted in the z-average diameter for this sample increasing to 63nm.

When one of the intensity size distributions shown in figure 1 is overlaid with the one shown in figure 2, the effect of the incorporation of the drug on the particle size of the microemulsion droplet can be clearly seen (figure 3). The peak due to the microemulsion droplet has increased in size due to the solubilisation of the drug molecules.

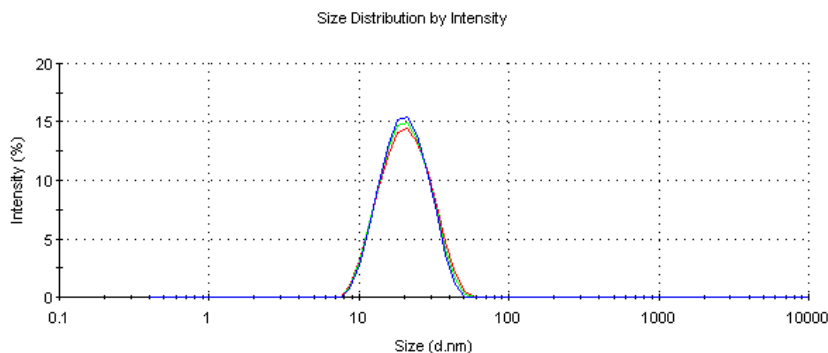
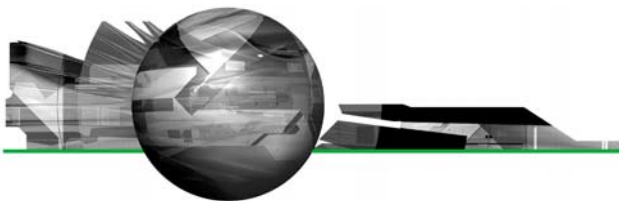


Figure 1: Intensity size distributions obtained from a microemulsion sample prepared by dilution of a pre-concentrate formulation in deionised water



Monitoring of the z-average diameter as a function of elapsed time since the formation of the microemulsion by dilution of the preconcentrate, allows the study of any loss of drug solubility. This information is invaluable in adjusting the formulation of the preconcentrate to maximize the solubility of the drug in the resulting microemulsion. The z-average diameter is intensity weighted and is therefore sensitive to the presence of large particles. It is a suitable parameter for following processes such as particle aggregation or crystallization.

Figure 4 shows the increase in the z-average diameter as a function of the elapsed time since the formation of the microemulsion by dilution of a preconcentrate containing an oil-soluble drug. This increase in the mean diameter results from the gradual loss of drug solubility from the microemulsion droplets. Suitability of various preconcentrates for maximizing drug solubility can be investigated by monitoring changes in the z-average diameter.

Conclusions

Dynamic light scattering is a non-invasive technique well suited to the study of colloidal dispersions such as microemulsions. The suitability of microemulsions for the solubilisation of drugs can be determined by following changes in the size distributions and mean diameters with time.

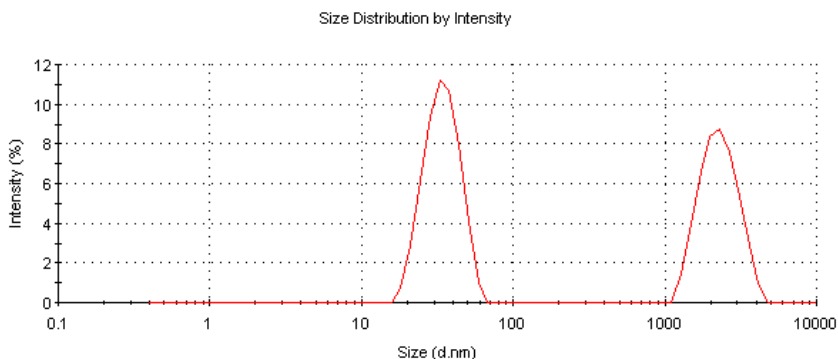


Figure 2: Intensity size distribution of a microemulsion obtained by dilution of a preconcentrate formulation containing an oil-soluble drug

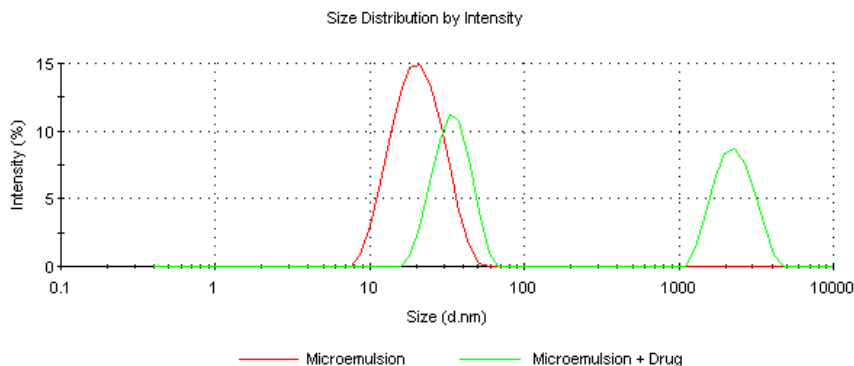


Figure 3: Overplot of the "empty" microemulsion and the microemulsion containing the oil-soluble drug

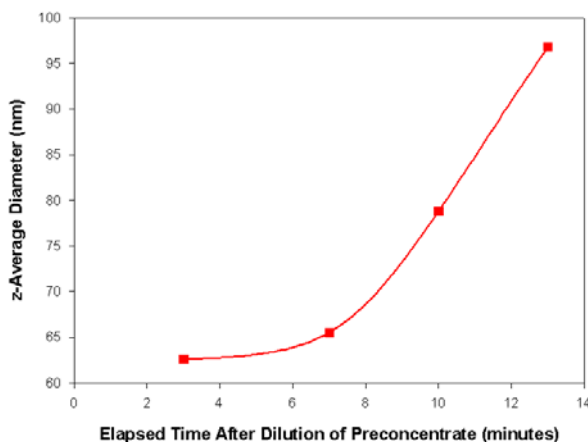
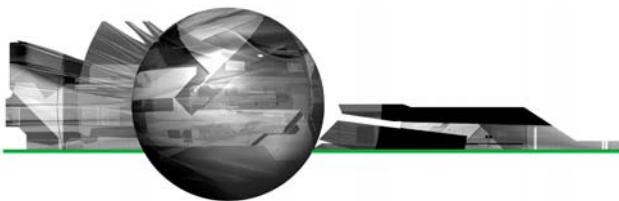


Figure 4: A graph showing the z-average diameter as a function of elapsed time since the formation of the microemulsion by dilution of the preconcentrate



Zetasizer Nano

The Zetasizer Nano system from Malvern Instruments is the first commercial instrument to include the hardware and software for combined static, dynamic, and electrophoretic light scattering measurements. The wide range of sample properties available for measurement with the Nano system include, particle size, molecular weight, and zeta potential.

The Zetasizer Nano system was specifically designed to meet the low concentration and sample volume requirements typically associated with pharmaceutical and biomolecular applications, along with the high concentration requirements for colloidal applications. Satisfying this unique mix of requirements was accomplished via the integration of a backscatter optical system and the design of a novel cell chamber. As a consequence of these features, the Zetasizer Nano specifications for sample size and concentration exceed those for any other commercially available dynamic light scattering instrument, with a size range of 0.6 nm to 6 μm , and a concentration range of 0.1 ppm to 40% w/v.

Complementing the patented hardware design, is the Malvern DTS software, providing instrument control and data analysis for the Zetasizer Nano System. The DTS software utilizes self analyzing algorithms to insure that the optical setup is optimized for each set of experimental conditions, and includes a unique "one click" measure, analyze, and report feature designed to minimize the new user learning curve.

Malvern Instruments Ltd

Enigma Business Park • Grovewood Road • Malvern • Worcestershire • UK • WR14 1XZ
Tel: +44 (0)1684 892456 • Fax: +44 (0)1684 892789

Malvern Instruments Worldwide

Sales and service centers in over 50 countries for details visit www.malvern.co.uk/contact

more information at www.malvern.co.uk