

## Instructions for Cat. #398 “NanoFuel-V Solvent” Diluent Solution

**NanoFuel-V Solvent** is NanoLight’s solution to the problem associated with the solubility of V-Coelenterazine and analogs thereof that have been made available recently. It differs from NanoLight’s original NanoFuel Solvent with its unique ability to dissolve very hydrophobic Coelenterazine analogues (like v-CTZ, MeO-v-CTZ-OMe or CTZ400a).

It is not designed for *IN VIVO* live animal imaging use; please use “Fuel-Inject-V” and sterile V-Coelenterazine for live animal imaging.



### Reconstituting V-Coelenterazine or its analogues using NanoFuel-V Solvent:

1. Warm NanoFuel-V to room temperature. (Use a dry heat block to speed up the process. Do not use a water bath. The liquid will draw moisture causing rapid degradation of v-Coelenterazine.)
2. Depending on the desired concentration of your stock solution add 500  $\mu$ l of NanoFuel-V Solvent to get a 1 mg/ml solution. Concentrations up to 4 mg/ml are possible.
3. Higher concentrations may need additional heating (60-80°C) for 5 min. Try to avoid using wet heat (see above).
4. This “stock-solution” is then added to buffer of choice for use in your experiments with a final concentration between 10 and 100  $\mu$ M.
5. Store the stock solution at -80°C. Avoid multiple freeze-thaw cycles. Minimize the time at elevated temperatures.
6. Do not use the solution if it turns dark brown/black. This is an indication of degradation (oxidation).