

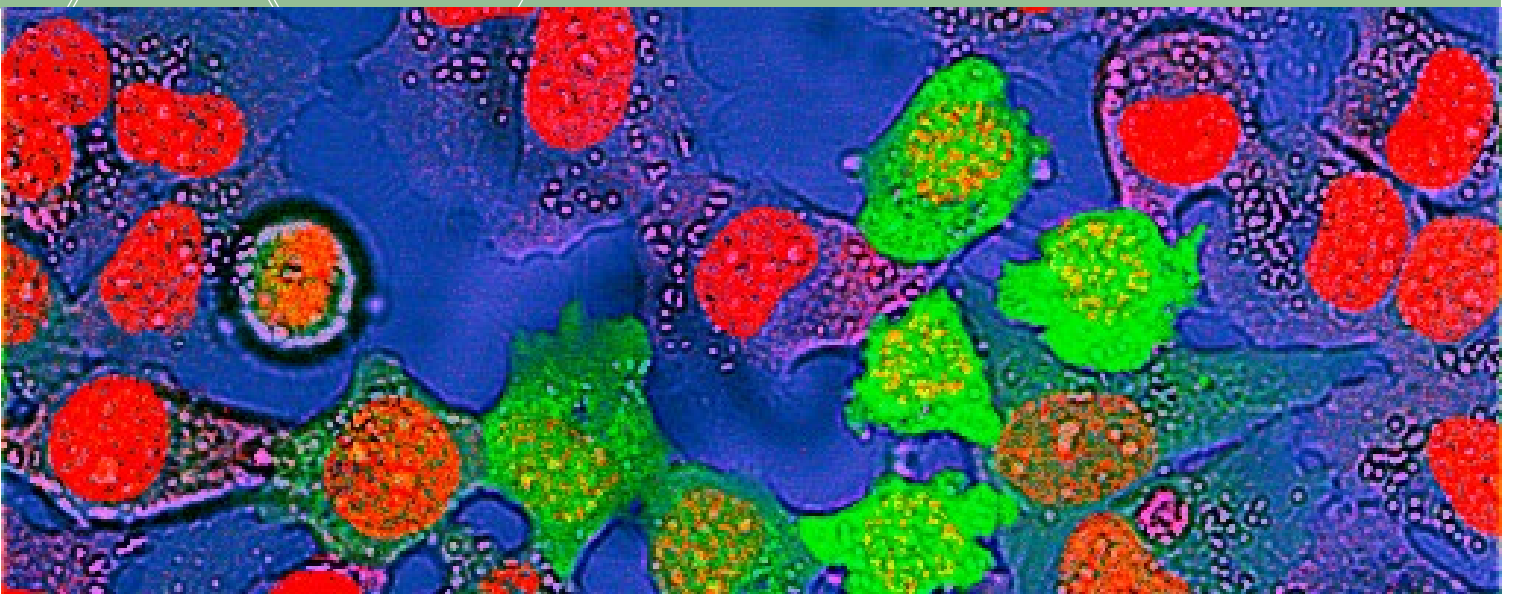
CyGEL™ SUSTAIN

A Novel Thermoreversible Mountant for Live Cells

IMAGING

CyGEL™

Analyse for LONGER with CyGEL SUSTAIN



- Specially formulated with a transition temperature for sol-to-gel at 24 degC enabling it to be handled as a sol at ambient temperature and a gel at 37 degC
- Formulated to allow the addition of RPMI and growth medium supplements for optimal analysis time - up to 2 hrs depending on cell type
- A medium formulation calculator is provided to allow users to establish customised culture media preparations for LIVE CELL preparations

CyGEL - it's a liquid when cold, and a gel when warmed

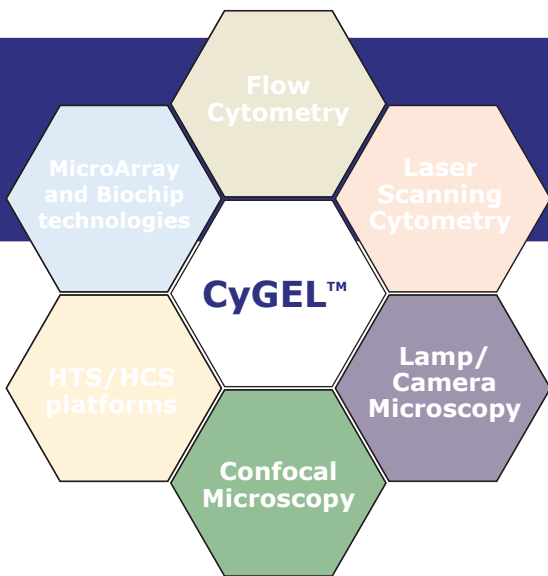
CyGEL SUSTAIN is supplied as a kit including:-

- > 10 x 500ul CyGEL SUSTAIN (specially formulated for sustained LIVE CELL analysis)
- > 1 x protocol
- > 1 x medium formulation calculator


Key Advantages:

- > A novel hydrogel compatible with LIVE Cells
- > Controllable and rapid transition from liquid to gel
- > Convenient immobilisation of LIVE Cells
- > Compatible with GFP and 'in gel' fluorescent probes including DRAQ5™
- > Optically clear with low autofluorescence

For more details please contact us, quoting product code:- CS20500



CyGEL™ SUSTAIN – a novel thermoreversible mountant for longer analysis of LIVE CELLS

CyGEL Sustain™ is manufactured by BioStatus Limited.  and distributed by ImmunoChemistry Technologies, LLC. Please contact:

ImmunoChemistry Technologies, LLC.
Toll Free: 800-829-3194



T: 952-948-0915 F: 952-888-8988
E: sales@immunochemistry.com W: www.immunochemistry.com

CyGEL™: MATERIAL SAFETY DATA SHEET. For further information see our website at www.biostatus.com

Ref: SO/230704/BS/CY15 **1. NUMBER(S)** CyGEL™ **2. COMPANY/UNDERTAKING NAME & ADDRESS** Biostatus Limited, 56 Charnwood Road, Shepshed, Leicestershire, LE12 9NP, United Kingdom **t:** +44 1509 558163 **f:** +44 1509 651061 **e:** enquiry@biostatus.com **w:** http://www.biostatus.com **3. COMPOSITION** Substance: CyGEL™. Chemical Description: POLOXAMER 407. Chemical Name: OXIRANE, METHYL-, POLYMER: CAS Number 9003-11-6 **4. HAZARDS IDENTIFICATION** This product is hazardous or contains components which are hazardous according to the OSHA Hazard Communication Standard. THIS PRODUCT MAY CONTAIN TRACE AMOUNTS OF ETHYLENE OXIDE, PROPYLENE OXIDE AND 1,4-DIOXANE, CHEMICALS KNOWN TO THE STATE OF CALIFORNIA TO CAUSE CANCER AND/OR BIRTH DEFECTS OR REPRODUCTIVE HARM. **5. FIRST AID MEASURES** Skin contact: WASH WITH SOAP & WATER; THEN FLUSH WITH WATER UNTIL CHEMICAL IS REMOVED. Remove and wash contaminated clothing before reuse. If irritation develops, get medical attention. Eye contact: EYE-FLUSH PROMPTLY WITH PLENTY OF COLD WATER FOR AT LEAST 15 MINUTES. If irritation develops, get medical attention. Ingestion: IF CONSCIOUS, INDUCE VOMITING AS DIRECTED BY DOCTOR, IMMEDIATELY GET MEDICAL HELP. Inhalation: REMOVE TO FRESH AIR. IF BREATHING IS DIFFICULT, GIVE OXYGEN. GET MEDICAL ATTENTION IF SYMPTOMS PERSIST. **6. ACCIDENTAL RELEASE MEASURES** Neutralizing Agent: COLD WATER. Methods for cleaning: WEAR PROTECTIVE EQUIPMENT. WIPE OR MOP UP AND PLACE IN A CONTAINER FOR DISPOSAL. FLUSH RESIDUE AWAY WITH COLD WATER. NON POROUS SURFACE CONTAMINATION MAY BE REMOVED USING COLD WATER **7. HANDLING AND STORAGE** Handling: LABORATORY COAT, RUBBER GLOVES AND SAFETY EYE GOGGLES Storage: STORE AT 4°C. DO NOT FREEZE. PROTECT CONTAINER FROM PHYSICAL DAMAGE. **8. EXPOSURE CONTROLS** Personal Protection Equipment: LABORATORY COAT. Eye protection: SAFETY GOGGLES. Hand protection: RUBBER GLOVES. **9. PHYSICAL AND CHEMICAL PROPERTIES** Appearance: CLEAR, TRANSPARENT SOLUTION. **10. STABILITY AND REACTIVITY** STABLE SOLUTION. NOT CORROSIVE, NON-OXIDISING, NO KNOWN HAZARDOUS DECOMPOSITION PRODUCTS. **11. TOXICOLOGICAL INFORMATION** Rabbit, Primary Skin Irritation: SLIGHTLY IRRITATING. Rat, Oral LD50: > 15g/Kg. PRACTICALLY NON-TOXIC. Rabbit, Corneal Anaesthetic Activity: NEGATIVE. NO CORNEAL ANAESTHETIC ACTIVITY. Rabbit, Assay for Pyrogenic Substances: NEGATIVE. PYROGEN FREE. Ames Salmonella Assay (Direct Plate): NEGATIVE, (Plate with S-9) NEGATIVE. NOT MUTAGENIC IN 5 STRAINS TESTED. Rabbit, Dermal LD50: > 15g/Kg: PRACTICALLY NON-TOXIC. Dog, Absorption through Open Wounds: NO ACCESS TO SYSTEMIC CIRCULATION. Mouse, intravenous LD50: 2.25g/Kg. SLIGHTLY TOXIC/PRACTICALLY NON-TOXIC. Rabbit, Eye Irritation (unrinsed) - (26% GEL): NON-IRRITATING. Acute Overexposure Effects: CONTACT WITH THE EYES AND SKIN MAY RESULT IN IRRITATION. INHALATION MAY RESULT IN RESPIRATORY IRRITATION. INGESTION MAY RESULT IN GASTRIC DISTURBANCES. Chronic Overexposure Effects: THERE ARE NO KNOWN CHRONIC EFFECTS ASSOCIATED WITH THIS MATERIAL. **12. DISPOSAL AND ENVIRONMENTAL CONSIDERATIONS** Spills or unwanted material should be placed in suitable containers for disposal in a licensed facility. Disposal should follow federal and local Regulatory Guidelines. **13. SOURCES OF KEY DATA** 1. BASF CORPORATION PERFORMANCE CHEMICALS <https://worldaccount.basf.com/wa/PublicMSDS> 2. BIOSTATUS LTD: <http://www.biostatus.co.uk> **14. LABEL DATA** Label Required: YES (TOXIC MATERIAL). Common Name: CyGEL™ Signal Word: WARNING! Acute Health Hazard: Moderate Contact Hazard: Slight Fire Hazard: Slight Reactivity Hazard: None Special Hazard Precautions: This product is hazardous or contains components which are hazardous according to the OSHA Hazard Communication Standard. THIS PRODUCT MAY CONTAIN TRACE AMOUNTS OF ETHYLENE OXIDE, PROPYLENE OXIDE AND 1,4-DIOXANE, CHEMICALS KNOWN TO THE STATE OF CALIFORNIA TO CAUSE CANCER AND/OR BIRTH DEFECTS OR REPRODUCTIVE HARM. Storage: STORE AT 4°C. DO NOT FREEZE. PROTECT FROM PHYSICAL DAMAGE.