

Liquid Scintillation Counters

Tri-Carb 4810TR

Low Activity Liquid Scintillation Counter



Tri-Carb 4810TR

Description

The Tri-Carb® 4810TR LSC is a computer controlled bench top liquid scintillation counter for detecting small amounts of alpha, beta and gamma radioactivity.

Standard instrument features

- Powerful built in computer system with Windows 8.1 operating system 4 GB (minimum) RAM and 250 GB hard disk (minimum) 3 high speed USB ports and Dual Gigabit Ethernet support with 19 inch Wide LCD monitor with DVI-D video output and keyboard.
- Date and time clock provides real time display and time-stamped printouts; battery supported.
- **Operational status indicator LED's.***
- Live SpectraView™ Automatic spectrum display aids in optimizing counting regions and helps evaluate complex sample situations. It allows setting temporary regions on the spectrum screen and enables the operator to monitor the effect of AEC (Automatic Efficiency Control) while the sample is counting.
- Quick-Count sample loading for 15 independent protocols (with unlimited assays) provides unrestricted access to sample changer and protocol selection plugs. Sample batches are processed by simply activating the Quick-Count protocol plugs, thus minimizing any user programming.
- Anti-jam recovery of the sample changer mechanism protects samples, vials and the counting system from damage if obstructions occur.
- Automatic power-fail recovery restarts counting when power is restored and the instrument has reinitialized itself.
- Positive sample identification provides protocol number, cassette number, sample number, and user-selectable printout and data file storage for the counting time and date on each sample.
- **A cassette-loaded bi-directional sample conveyor mechanism is standard with a sample capacity of either 408 standard 20 mL vials, or 720 small 4 or 7 mL vials.***
- Multi-parameter linear MCA (Multichannel Analyzer) with an effective resolution of 1/10 keV, offers an extended dynamic quench range and provides multi-parameter spectrum analysis to correct for luminescence, color quenching and background radiation.
- **Patented TR-LSC® (Time-Resolved Liquid Scintillation Counting) is featured for high sensitivity, low background counting of LS samples. TR-LSC increases sample throughput.***

- ^{133}Ba low energy external standard source and tSIE (transformed Spectral Index of External standard). The use of integral spectrum counts eliminates the need for repeat counting of the external standard and negates the effect of isotope on quench monitoring accuracy and precision. The ^{133}Ba external standard is centered under the sample vial which eliminates the effects of volume variations and assures reproducible quench monitoring for the life of the instrument.*
- User adjustable assay-specific sample precount delay permits dark adaptation of samples before counting.

Standard software features

- QuantaSmart™ software with comprehensive on-line context sensitive help for the Windows® 8 operating system is a 32-bit operating system software that provides a robust multitasking, easy networking environment and unlimited assays in a secure multiuser environment.*
- Enhanced Direct DPM technique determines the DPM of any single label pure beta or beta/gamma radionuclide in any cocktail without the use of quench standards.
- Luminescence detection and reporting with percent luminescence is flagged on printout to alert user of possible sample problems.
- Assay-specific, user selectable, coincidence resolving time enables optimized counting of any fast or slow, liquid or solid scintillator, for SPAs (Scintillation Proximity Assays), solid scintillation filters and the newer liquid scintillation cocktails.
- Spectral unfolding separates and displays the individual radionuclide spectra of dual label samples in color (requires color-corrected dual label DPM option) analysis of sample spectrum.
- 3D (three-dimensional) spectral mapping displays in color the quench standard spectra together with the spectrum of the unknown for single label DPM counting (requires color-corrected dual label DPM option).
- SpectraBase counting and data management system provides region less counting and storing of complete spectra for all samples and standards. Features include automatic recall of spectra stored in the quench library for region-independent quench correction and post processing of sample data with the optional Replay™ feature.*
- Decay computations automatically calculate decay corrected DPM values for commonly used radionuclide standards.
- Group PrioStat™ interrupt mode gives priority counting status to a batch of samples counted according to any stored protocol conditions. It automatically restores the interrupted protocol upon completion and stores PrioStat data for immediate viewing. Data is printed at protocol termination.

- Background subtraction can be nominated via sample, entered value or stored IPA background spectrum.
- SIS (Spectral Index of Sample) quench indicating parameter determines counting efficiency by analysis of sample spectrum.
- Programmable single photon counting enables luminescence assay counting with optimized signal-to-background ratios. It overcomes problems associated with excessive luminescence.
- Preset time (up to 9,999.99 minutes) and preset error coincidence termination optimizes counting accuracy in the three counting regions.
- Automatic spectrum plot (on demand) per sample allows spectral documentation of samples.
- Sample screening allows screening numeric fields such as activity based on several criteria including background levels, a hard number or within a range of activities or values. Hits can easily be identified in reports with optional highlighting and custom hit flags.
- Printed header contains instrument serial number, user ID, and drive and path of all electronic stored data. Each printed page or RTF (Rich Text Format) file report is numbered and dated for GLP compliance.
- Password protection prevents unwanted changes to saved assays.
- Half-life correction to any date and time is available for up to three radionuclides.
- Activity reporting is provided in becquerels, micro-curies, or picocuries.
- Auto QA (Automatic Quality Assurance) automatically prints reports for backgrounds, efficiencies, E2/B, and Chi-square values for both ^3H and ^{14}C including the date and time for each test. Auto QA results may be transmitted via RS-232 for archiving.
- Percent of standard calculations is present for single, dual and triple label samples.
- Automatic processing of count data to final results provides automatic, protocol specific data processing for all user applications, including commercial or user-generated software. No exporting of data to off board storage devices or computers is required.
- Independent output formatting to printer, for each protocol provides almost unlimited flexibility in data reporting. Electronic data can be saved to disk in ASCII, RTF, or Microsoft® Excel® compatible formats. Reports can be customized for data content and protocol information.
- Computer-aided diagnostics are used to verify all system functions.

Hardware options

- Automatic 2D barcode reader reads 2D barcodes (factory default setup). Barcode readings can be used to create the sample work list (work list option required) and optionally save to a file or validate an existing work list. Barcodes are enabled on individual assays giving the user maximum flexibility in barcode usage.*
- Varisette™ sample changer enables intermixing and counting of both large and small sample vials without special adapters. Option includes both large vial (12-position) and small vial (18-position) cassettes.*
- Printer (ink jet or laser jet).
- Temperature-controlled refrigeration establishes and maintains optimum counting conditions for a wide variety of sample types.

Software options

- Sample work list enables entry, editing and review of work lists for each assay. Automatic creation of work lists is possible with the 2D barcode option, which allows sample identification with user-specified codes for sample printouts and data files.
- Dynamic color-corrected single and dual label DPM is based on tSIE with AEC. AEC corrects for differential quenching effects in multi-label samples. The low energy spectrum of the external standard ensures accurate tracking of ^3H , ^{14}C and other low energy sample spectra over a very wide quench range. Includes factory stored quench standards for ^3H and ^{14}C in classical and ULTIMA Gold™ cocktails with On-screen editing of quench correction curves and recording of date last modified.*
- Enhanced IPA™ (Instrument Performance Assessment) database for monitoring efficiencies, backgrounds, E2/B and Chi-square values for ^3H and ^{14}C over the life of the instrument. IPA flags impending problems and provides both running mean and fixed baseline charts and associated tables for retrospective quality control and proactive system maintenance. Baseline acquisition is programmable for increased flexibility. The most recent IPA time and date stamped data are available on demand for reporting purposes. Each IPA printout includes instrument model, serial number, software version number and calibration standard information.*
- Enhanced Replay™ sample post-processing provides complete recall and post-processing of historical count data to eliminate sample recounting. It enables changes to count conditions and reports as well as execution of user application software for optimization of data analysis.*

- HSCM (High Sensitivity Count Mode) increases system sensitivity by implementing additional electronic background reduction via TR-LSC. It includes assay specific, user-selectable delay before burst settings for optimization of TR-LSC.
- Triple-label DPM is based on tSIE/AEC for accurate spill correction (requires color-corrected single and dual label DPM option).
- Chemiluminescence correction with response normalization corrects for luminescence interference to speed up sample counting. Response normalization of the correction circuits eliminates the effect of component drift on the corrected results.
- SpectraWorks™ spectrum analysis software for the Windows® operating system analyzes beta, alpha, and gamma spectra and provides simultaneous display for up to four spectra in stacked or overlaid mode. It features zooming to any part of the spectrum; six regions of interest; display of counts or CPM and linear or log spectra; provides automatic and manual scaling; calculates E2/B, MDA, peak resolution; allows adding and subtracting of spectra and multiplication and division by constants.
- Easy View Raw data saved on the PC hard disk or a network drive can later be processed off-line with EASY View spectrum analysis software. EASY View displays up to six spectra simultaneously and allows spectral arithmetic, DPM calculations, statistical analyses and radiocarbon age dating.
- 60-user protocol option upgrades the 4810TR to allow for 60-user programming capability.

Accessories

- Instrument utility cart: functionally designed general purpose laboratory cart. Supports any PerkinElmer bench top system.
- See the Equipment, Chemicals & Supplies section in the PerkinElmer catalog.

Physical data

Dimensions:

Height: 18.5 in. (47 cm)

Width: 40.5 in. (103 cm)

Depth: 32 in. (81 cm)

Depth with refrigeration: 44 in. (112 cm)

Weight:

477 lb (217 kg) with refrigeration

Shipping weight: approximately 700 lb (318 kg)

Electrical Requirements:

100-240 Vac 50/60 Hz

3-prong grounded plug

Power Consumption:

<200 VA; <800 VA with temperature control option

Environmental:

Operating ambient temperature 15 to 35 °C (59 to 90 °F)

Operating relative humidity 30% to 85%

Typical performance data

(As measured at Singapore facility.)

Energy Range: 0–2,000 keV

Efficiency, Normal Count Mode:

| | | Minimum Acceptable |
|-----------------|------------|--------------------|
| ³ H | 0–18.6 keV | 60% |
| ¹⁴ C | 0–156 keV | 95% |

Figure of Merit (E²/B), Normal Count Mode (NCM):

| | | |
|-----------------|------------|-----|
| ³ H | 1–18.6 keV | 180 |
| ¹⁴ C | 4–156 keV | 380 |

Figure of Merit (E²/B), Low Activity/High Sensitivity Count Mode (HSCM):

| | | |
|-----------------|---------------|-----|
| ³ H | 1–12.5 keV | 300 |
| ¹⁴ C | 14.5–97.5 keV | 950 |

Observed Background, NCM:

| | | Average |
|-----------------|------------|---------|
| ³ H | 0–18.6 keV | 17 CPM |
| ¹⁴ C | 0–156 keV | 26 CPM |

Note: The efficiencies, backgrounds, and E²/B values for the Normal Count Mode were determined using using PerkinElmer sealed large vial glass standards set P.N. 6008500 verified with NIST standard activity.

The HSCM values are determined using PerkinElmer low level sealed large glass vial standards set P.N. 6008914 verified with NIST standard activity. No maximum is specified for background. Safety, Radiated Emissions and Immunity:

The Tri-Carb 4810TR has been tested and approved for safety, radiated emissions and immunity according to the standards of UL, IEC61010 and CE.

In the U.S.A. the UL approval satisfies the requirements of 29CFR 1910.399.

*Items marked with an asterisk are unique and exclusive features or specifications to the PerkinElmer Liquid Scintillation Instruments.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2009-2015, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

007089E_01

PKI