



## PERKINELMER INFORMATICS ADVANCED CONNECTOR FOR SAS®

### Faster Answers with Smarter Workflows

Our Advanced Connector for SAS® provides a high performance and scalable solution over the current out-of-the-box integration between TIBCO Spotfire® and SAS®. The connector is built on a modern web service framework that is faster and more scalable compared to traditional OLE DB provider based loads. The Advanced Connector for SAS® empowers users to create more robust applications, data flows and integrations with TIBCO Spotfire® that have dependencies on SAS® source datasets and transport files.

### PerkinElmer Informatics Advanced Connector for SAS®



PerkinElmer is the exclusive global distributor of the TIBCO Spotfire® platform for certain scientific and clinical R&D applications.

### AT A GLANCE

- Connectivity with SAS Datasets and Transport files
- High-Performance connectivity via SAS BI Web Services
- Out-of-the-box CDISC SDTM Knowledge
- Effortlessly pool multiple study-data into one analysis from multiple sources
- Self-documenting data transactions
- Low-Maintenance Security (role-based account privilege assignment)
- Intuitive User Experience

**High Performance, Scalability, and Out-of-the-Box Knowledge of CDISC SDTM delivered via Web Services based Framework**

### RETURN ON INVESTMENT FROM CUSTOMERS:

- Faster review of clinical data across all studies, accelerating time to market for products across the entire portfolio of clinical candidates.
- Ability to build robust and enterprise-scale applications using a combination of TIBCO Spotfire® and SAS datasets
- Ensure Data Standard compliance while providing an easy way to adhere to data standards
- Lower IT involvement in data access assignment and maintenance of data flows

# ATTRIBUTES AND CAPABILITIES

## Out-of-the-box CDISC SDTM Knowledge

The screenshot displays the 'Clinical/Translational Administration' software interface. On the left, a tree view shows the 'SDTM Studies - Datasets' structure, including 'Program001\CDISC PILOT01' and 'Program003\study01'. The 'AE - Adverse Events [AE]' dataset is selected. The main window shows the 'Specification Name' and 'Spottier Column Name' mapping table for 'AE - Adverse Events [AE]'. The table lists various AE-related terms and their corresponding data types and connector types.

Specification Name	Spottier Column Name	Data Type	Source Field	Connector Type	Matched By
AEACN	Action Taken with Study Treatment	Text	AEACN	Text	Auto
AEACNTH	Other Action Taken	Text			
AEBDSYCD	Body System or Organ Class Code	Numeric			
AEBDSYS	Body System or Organ Class	Text	AEBDSYS	Text	Auto
AECAT	Category for Adverse Event	Text			
AECTRTR	Concomitant or Additional Treatment Given	Text			
AEDECOD	Dictionary-Derived Term	Text	AEDECOD	Text	Auto
AEDUR	Duration of Adverse Event	Text			
AEENDTC	End Date/Time of Adverse Event	Date/Time	AEENDTC	Text	Auto
AEENDY	Study Day of End of Adverse Event	Integer	AEENDY	Numeric	Auto
AEENRF	End Relative to Reference Period	Text			
AEENRPT	End Relative to Reference Time Point	Text			
AEENPT	End Reference Time Point	Text			
AERPID	Group ID	Text			
AEHLGT	High Level Group Term	Text			
AEHLGTC	High Level Group Term Code	Numeric			
AEHLT	High Level Term	Text			
AEHLTCD	High Level Term Code	Numeric			
AELLT	Lowest Level Term	Text			
AELLTCD	Lowest Level Term Code	Numeric			
AELOC	Location of Event	Text			
AEMODIFY	Modified Report Term	Text			

## Advanced Connector for SAS® populating PerkinElmer's Solution for Clinical Data Review

The screenshot displays the 'Advanced Connector for SAS' interface. The top section shows a horizontal bar chart titled 'Adverse Events - Patient Count by Study Arm'. The chart lists various adverse event categories and their corresponding patient counts. The categories are color-coded by severity: MILD (green), MODERATE (yellow), and SEVERE (red).

Adverse Event Category	Count
GENERAL DISORDERS AND ADMINISTRATION SITE CONDITIONS	108
SKIN AND SUBCUTANEOUS TISSUE DISORDERS	106
NERVOUS SYSTEM DISORDERS	59
GASTROINTESTINAL DISORDERS	53
CARDIAC DISORDERS	44
INFECTIONS AND INFESTATIONS	39
RESPIRATORY, THORACIC AND MEDIASTINAL DISORDERS	30
PSYCHIATRIC DISORDERS	29
INVESTIGATIONS	23
MUSCULOSKELETAL AND CONNECTIVE TISSUE DISORDERS	20
INJURY, POISONING AND PROCEDURAL COMPLICATIONS	14
RENAL AND URINARY DISORDERS	11
METABOLISM AND NUTRITION DISORDERS	10
VASCULAR DISORDERS	8
EYE DISORDERS	7
SURGICAL AND MEDICAL PROCEDURES	5
EAR AND LABYRINTH DISORDERS	4
CONGENITAL, FAMILIAL AND GENETIC DISORDERS	3
NEOPLASMS BENIGN, MALIGNANT AND UNSPECIFIED (INCL. CYSTS AND...)	3
REPRODUCTIVE SYSTEM AND BREAST DISORDERS	3
IMMUNE SYSTEM DISORDERS	2
HEPATOBIILIARY DISORDERS	1
SOCIAL CIRCUMSTANCES	1

The bottom section shows a heatmap titled 'Adverse Events - Patient Count by Study Arm'. The heatmap displays the distribution of adverse events across different study arms (Xan\_Hi, Xan\_Lo, Pbo) for various body systems and organ classes. The cells are color-coded based on the patient count, with green representing low counts and red representing high counts.

To learn more about our solutions for clinical development, please visit [www.perkinelmer.com/ClinicalInformatics](http://www.perkinelmer.com/ClinicalInformatics)

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

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

012562\_01

PKI