

I soci devono documentare e rendere pubblici ogni anno attraverso un apposito Modello di Trasparenza che costituisce parte integrante del presente Codice, i trasferimenti di valore effettuati direttamente o indirettamente ai Professionisti del Settore Sanitario, alle Organizzazioni Sanitarie e alle Terze Parti.

	Full Name	HCPs: City of Principal Practice HCOs: City where registered	Country of Principal Practice	Principal Practice Address	Donations	Scholarship	Contribution to costs of Events			Contribution to costs of Training			Costs for the participation in training, educational and promotional events on company's products organised by the Members	Fee for service and consultancy		Transfers of Value Research & Development	Total	
							Sponsorship agreements with HCOs/Third Parties appointed by HCOs to manage an event	Registration Fees	Travel & Accommodation	Sponsorship agreements with HCOs/Third Parties appointed by HCOs	Registration Fees	Travel & Accommodation		Fees	Related expenses agreed in the fee for service or consultancy contract, including			
HCPs (Healthcare Professionals)	INDIVIDUAL NAMED DISCLOSURE - one line per HCP (i.e. all transfers of value during a year for an individual HCP will be summed up: Remitization should be available for the individual Recipient or public authorities' consultation only, as appropriate)																	
	Daniela Verginelli	Roma	Italy	Istituto Zooprofilattico Sperimentale Lazio e Toscana M. Aleandri - Via Appia Nuova,1411	NA	NA	NA*	NA*	NA*	NA	NA	NA	NA	1300	NA	1300		
	Pr. Vittorio Sambri	Bologna	Italy	Via Odone Belluzzi 9, 40135	NA	NA	NA	NA	NA	NA	NA	508,8	NA	NA	508,8			
	Dr. Luca Falzone	Catania	Italy	Piazza Università, 2, 95124 Catania CT	NA	NA	NA	NA	NA	NA	NA	NA	750,4	NA	750,4			
	Dr. Giovanna Travi	Segrate	Italy	Via Siena 2/4	NA	NA	NA*	NA*	NA*	NA	NA	NA	600	NA	600			
	AGGREGATE DISCLOSURE - per HCPs																	
	Aggregate amount attributable to transfers of value to such Recipients					NA	NA	NA*	NA*	NA*	NA	NA	NA	NA	800	800		
Healthcare Third Parties	INDIVIDUAL NAMED DISCLOSURE - one line per HCOs (i.e. all transfers of value during a year for an individual HCOs will be summed up: Itemization should be available for the individual Recipient or public authorities' consultation only, as appropriate)																	
	Kassiopea Group Srl	Cagliari CA	Italy	Torre Verde, Via S. Tommaso D'Aquino, 20, 09134	Only HCO	NA	1200	NA	NA	NA	NA	NA	NA	NA	1200			
	VET INTERNATIONAL SRL	MILANO	Italy	Via Carlo Farini 81	Only HCO	NA	4000	NA	NA	NA	NA	NA	NA	NA	4000			
	SIMTI PRO	MILANO	Italy	Via Desiderio, 21, 20131	Only HCO	NA	20740	NA	NA	39711	NA	NA	NA	NA	60451			
	SYMPOSIUM MEETING E CONGRESS	Catanzaro CZ	Italy	Via Raffaele Teti, 79, 88100	Only HCO	NA	3660	NA	NA	NA	NA	NA	NA	NA	3660			
	Meeting srl	Udine	Italy	Via Villalta 32	Only HCO	NA	9760	NA	NA	NA	NA	NA	NA	NA	9760			
	AFFIDIA SRL SB	Trieste	Italy	Località Basovizza 558	Only HCO	NA	3500	NA	NA	NA	NA	NA	NA	NA	3500			
Healthcare Organisations	AGGREGATE DISCLOSURE																	
	Aggregate amount attributable to transfers of value to such Recipients				Only Third Parties	NA	NA	NA	NA	NA	NA	NA	NA	NA				
	Number of Recipients in aggregate disclosure				Only Third Parties	NA	NA	NA	NA	NA	NA	NA	NA	NA				
	% of the number of Recipients included in the aggregate disclosure in the total number of Recipients disclosed				Only Third Parties	NA	NA	NA	NA	NA	NA	NA	NA	NA				

* In the case of direct support for the formation of HCPs that they exercise in a private context, it will be necessary to publish the data in individual or aggregate form, depending on whether the HCP has given consent